

life technologies™

Life Technologies
850 Lincoln Center
Foster City, CA 94404
Gregory.Govoni@lifestech.com

Applied Biosystems | **Invitrogen**

¹ Applied Biosystems, 850 Lincoln Centre Drive, Foster City, CA, USA; ² Cornell University, Ithaca, USA.

New generation sequencing technology provides researchers access to unprecedented amounts of genetic information. This genetic information can be parsed and analyzed for any number of disease determinants, including microbial pathogenicity. In this study, we sequenced whole genomes from 16 different serotypes of *Salmonella enterica* subsp. *enterica* representing 13 different serotypes (D2, E1, E4, F, G, I, L, N, O, O54, P, Q, and R). Mate-pair libraries for each genome were created and analyzed using the SOLiD™ sequencing system. From the data generated, a de novo assembly for each genome was constructed and analyzed for the presence of all 21 reported *Salmonella* Pathogenicity Islands (SPIs) as well as several LT2 prophages (fels1- & 2, gifsy-1 & 2). For many of the SPIs, their presence was universal (SPIs 1-6, 9, 11-13). SPIs that were not found in this subset of strains, and seem to be serotype-specific include SPIs 10, 15, 17, and 19-21. Frequency of the LT2 prophage insertions varied between 31% and 62% of the strains analyzed. Further analysis finds that only the fels-2 insertion has been maintained in a similar context as in the LT2 genome. The other remaining prophage insertions have weak identity scores and lack consistent genomic context, which suggests these insertions may have occurred independently with distantly-related prophages. We also took a more detailed view of the genes that make up the SPIs. First, we examined the presence or absence of several SPI genes that have been shown to be important for virulence. Interestingly, several orthologues for known virulence factors (*avjA*, *sseA*, *sseB* and *STM1318*) were missing in at least 1 isolate. Second, we looked at sequence variations in several of these SPI virulence factors orthologues. While some had elevated overall distance averages (Dn), *sipD* had the highest variation with a Dn of 0.0543 (as compared to 0.004-0.006 for highly conserved genes). Preliminary phylogenetic analysis grouped these sequences into several clusters. Interestingly for the *sseD* orthologues, the Uganda and Senftenberg isolates shared their own independent cluster, while the Uganda and Give isolates clustered separately despite being from the same E1 serogroup. Since the Uganda and Senftenberg isolates both came from a bovine source and the Give isolate did not (human), these findings may have possible implications for discerning the molecular basis behind *Salmonella* strain host specificity.

Salmonella enterica is a foodborne pathogen which causes an estimated 1.4 million human cases annually in the US [Mead et al., 1999]. It is a highly diverse pathogen with more than 2,500 different serotypes of *Salmonella* described. The host range and disease severity varies depending on the serovar in question. For example, *Salmonella enteritidis* causes a limited gastroenteritis in a broad range of hosts, while *Salmonella typhi* can cause a systemic, typhoid fever disease specifically in primates. Many of the genes responsible for these variations in host range and disease severity are due to virulence factors found in *Salmonella* Pathogenicity Islands (SPIs). SPIs are genomic regions in the chromosome that have been horizontally acquired and are important for disease pathogenesis in one form or another. To date, 21 SPIs have been described in the literature. While many islands have been studied in-depth (such as SPI-1 and SPI-2), many SPIs have only been recently reported and are not well-characterized. In general, not much is known regarding the conservation of these SPIs across a broad range of serogroups. The development of new molecular-based technologies is making a comprehensive analysis of SPIs easier and faster to do. SOLID™ next generation sequencing is an incredibly powerful tool that allows researchers to analyze a vast amount of genomic data in a relatively quick amount of time. Now it is possible to look in parallel at every SPI in a diverse set of strains in a matter of months without sacrificing on the level of detail required to do real in-depth analysis. In the study reported here, we describe the prevalence of each of these SPIs in the strains tested and perform some of the analyses that can be done to help discern which genes may play a role in determining host specificity and/or disease severity.

Table 1. List of Strains Used in this Study

Source	Host	Serogroup	Source	Strain
100	human	O:6:H:3	97-6	Atlanta
101	human	O:6:H:3	97-13	Atlanta
102	human	O:4:H:3	98-1	Atlanta
103	human	O:16:H:1	99-17	Genitum
104	human	O:13:H:1	99-15, 99-16, 99-17	Genitum
105	human	O:15:H:1	99-18	HaitianUS
106	human	O:3:H:1	99-19	horne
107	vine	O:15:H:1	1-2000-01	horne
108	human	O:2:H:1	21-00-01	Mexico
109	human	O:2:H:1	21-00-02	Mexico

Genome sequencing and assembly. Genomes were sequenced using the SOLiD™ system (Applied Biosystems, Foster City). Mate-paired libraries with approximately 1.5 kb inserts were constructed and deposited on one-eighth of a flowcell. Then, 25 bp reads were obtained from each of the F3 and R3 tags. After correcting errors in colospace reads using the modified version of the spectral alignment tool called the EULER-USR package (Chaisson, et al., 2009), *de novo* assembly was performed using the SOLiD™ *de novo* pipeline, which employs the Velvet assembly engine (Zerbino & Birney, 2006).

Scaffolds were aligned to two reference genomes (*S. Typhimurium* LT2 and *S. Enteritidis*) and concatenated into pseudogenomes. Scaffolds that did not match the chromosomes of the reference genomes were considered to be putative plasmids or strain specific transposable elements. Read coverage was approximately 120–180 fold per nucleotide.

Whole genome and SPI Sequence Analysis. Automated annotation was performed with the RAST server (<http://rast.nmpdr.org>; Aziz et al. 2008). The reference sequences used to delineate the SPIs and LT2 prophages are listed in Table II and were obtained from multiple sources such as the PAI database (Yoon et al. 2007). Individual ORF analysis was performed by taking ORF sequences from reference strain and generated the *Salmonella* LT2 reference assemblies with BLASTN. To ensure basic hits were in the proper genomic context, the sequence based comparison tool on the RAST server was used to align the pseudogenomes against selected reference genomes.

References	glyA2 prophage	S. typhimurium (B2)	45.8	48	1080228	STM1005	STM1052	1145450
Aziz RK, Bartels D, et al. 2008. The RAST Server: Rapid Annotations using Subsystems Technology. BMC Genomics 9(7): 336-46.								
Chaisson MJ, Brinza D, Pevzner PA. 2009. De novo fragment assembly with short mate-paired reads: Does the read length matter? Genome Res 19(2): 336-46.								
Darling AC, Ma B, Blattner FR, Perna NT. 2004. De novo: multiple alignment of conserved genomic sequence with rearrangements. Genome Res 14(7):1394-401.								
Dickins JE, Jernett E. 2002. Velvet: algorithms for de novo short read assembly using the paired-end approach. Genome Res 12(12):1812-19.								
Mead, P. S., Lutskevich, V., Dietz, L. F., McCaig, J. S., Breesee, C., Shapiro, P. M., Griffin, and R. V. Tauxe. 1999. Food-related illness and death in the United States. Emerg. Infect. Dis. 5:607-625.								
Yoon SH, Park YK, Lee S, Cho D, HK TK, Cho K, Kim JF. 2007. Towards Pathogenomics: A Web-Based Resource for Pathology Analysis. Nucleic Acids Research 35 D395-D400.								

For Research Use Only. Not intended for any animal or human therapeutic or diagnostic use.
© 2010 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation and their respective owners.

[illegible]

Figure 1. Legend: A combination of blackies and the sequence based comparison tool on the RAST server was used to determine the presence or absence of every coding sequence (CDS) in every SPI. The % of CDSs present in each isolate relative to the number in the reference strain are reported for each SPI. Isolates with CDSs present in the SPI but do NOT appear to share the same genomic context as the reference strain are denoted with a box around the %. Strains with similar CDS order but with low CDS homology have percentages in purple. **Legend:** All the strains tested and include SPIs 1-6, 9, 11-13. Of those 11 SPIs, SPIs 1, 2, 4 and 5 are retained with over 90% of the reference ORFs present, suggesting there is evolutionary pressure to keep most of those islands intact. Islands that are not present in any of the isolates and can be classified as highly serovar-specific include SPIs 10, 15, and SPIs 19-21 are also most likely serovar specific, but may be present in these strains due to an independent horizontal acquisition of a related piece of DNA. Strains with SPIs 14 and 17 are the only two of the strains tested but located in different regions of the genome. Strains with significant portions of SPI7 include Inverness, Rubislaw, and Urbana.

A.

Vs. LT2

Ad - Outside
Ala
Ba
Cm
Gv
Hv
In
Jo - Inside

γ f-1
γ g-1

γ f-2
γ g-2

p

B.

Vs. CT18

γ 1
γ 2
γ 3

4 f
6 p

Vs. LT2

Min - Outside
Ms
Mon
Ru
Se
Ug
Urb
W - Inside

γ f-1
γ g-1

γ f-2
γ g-2

p

4 f
5 p
6 p

γ 1
γ 2
γ 3

▲ = potential prophage remnant

Percent protein sequence identity

Bidirectional best hit: 100 99.9 99.8 99.5 99 98 95 90 80 70 60 50 40 30 20 10

Unidirectional best hit: 100 99.9 99.8 99.5 99 98 95 90 80 70 60 50 40 30 20 10

Fig. 2. Legend: The sequence-based comparison tool on the RAST servers was used to generate color-coded comparative representations of the *S. Typhimurium* LT2 (A) or *S. typhi* CT18 (B) genome versus the pseudogenomes from the 16 newly sequenced isolates. Each putative coding sequence (CDS) is color-coded according to its percent protein identity to the reference genome. In Fig. 2A, the L2-4 prophage insertions and several other genes (e.g. *hlyD*, *hlyE*, *hlyF*, *hlyN*, *hlyP*, *hlyQ*, *hlyR*, *hlyS*, *hlyT*, *hlyU*, *hlyV*, *hlyW*, *hlyX*, *hlyY*, *hlyZ*, *hlyA*, *hlyB*, *hlyC*, *hlyD*, *hlyE*, *hlyF*, *hlyG*, *hlyH*, *hlyI*, *hlyJ*, *hlyK*, *hlyL*, *hlyM*, *hlyN*, *hlyO*, *hlyP*, *hlyQ*, *hlyR*, *hlyS*, *hlyT*, *hlyU*, *hlyV*, *hlyW*, *hlyX*, *hlyY*, *hlyZ*, *hlyA*, *hlyB*, *hlyC*, *hlyD*, *hlyE*, *hlyF*, *hlyG*, *hlyH*, *hlyI*, *hlyJ*, *hlyK*, *hlyL*, *hlyM*, *hlyN*, *hlyO*, *hlyP*, *hlyQ*, *hlyR*, *hlyS*, *hlyT*, *hlyU*, *hlyV*, *hlyW*, *hlyX*, *hlyY*, *hlyZ*, *hlyA*, *hlyB*, *hlyC*, *hlyD*, *hlyE*, *hlyF*, *hlyG*, *hlyH*, *hlyI*, *hlyJ*, *hlyK*, *hlyL*, *hlyM*, *hlyN*, *hlyO*, *hlyP*, *hlyQ*, *hlyR*, *hlyS*, *hlyT*, *hlyU*, *hlyV*, *hlyW*, *hlyX*, *hlyY*, *hlyZ*, *hlyA*, *hlyB*, *hlyC*, *hlyD*, *hlyE*, *hlyF*, *hlyG*, *hlyH*, *hlyI*, *hlyJ*, *hlyK*, *hlyL*, *hlyM*, *hlyN*, *hlyO*, *hlyP*, *hlyQ*, *hlyR*, *hlyS*, *hlyT*, *hlyU*, *hlyV*, *hlyW*, *hlyX*, *hlyY*, *hlyZ*, *hlyA*, *hlyB*, *hlyC*, *hlyD*, *hlyE*, *hlyF*, *hlyG*, *hlyH*, *hlyI*, *hlyJ*, *hlyK*, *hlyL*, *hlyM*, *hlyN*, *hlyO*, *hlyP*, *hlyQ*, *hlyR*, *hlyS*, *hlyT*, *hlyU*, *hlyV*, *hlyW*, *hlyX*, *hlyY*, *hlyZ*, *hlyA*, *hlyB*, *hlyC*, *hlyD*, *hlyE*, *hlyF*, *hlyG*, *hlyH*, *hlyI*, *hlyJ*, *hlyK*, *hlyL*, *hlyM*, *hlyN*, *hlyO*, *hlyP*, *hlyQ*, *hlyR*, *hlyS*, *hlyT*, *hlyU*, *hlyV*, *hlyW*, *hlyX*, *hlyY*, *hlyZ*, *hlyA*, *hlyB*, *hlyC*, *hlyD*, *hlyE*, *hlyF*, *hlyG*, *hlyH*, *hlyI*, *hlyJ*, *hlyK*, *hlyL*, *hlyM*, *hlyN*, *hlyO*, *hlyP*, *hlyQ*, *hlyR*, *hlyS*, *hlyT*, *hlyU*, *hlyV*, *hlyW*, *hlyX*, *hlyY*, *hlyZ*, *hlyA*, *hlyB*, *hlyC*, *hlyD*, *hlyE*, *hlyF*, *hlyG*, *hlyH*, *hlyI*, *hlyJ*, *hlyK*, *hlyL*, *hlyM*, *hlyN*, *hlyO*, *hlyP*, *hlyQ*, *hlyR*, *hlyS*, *hlyT*, *hlyU*, *hlyV*, *hlyW*, *hlyX*, *hlyY*, *hlyZ*, *hlyA*, *hlyB*, *hlyC*, *hlyD*, *hlyE*, *hlyF*, *hlyG*, *hlyH*, *hlyI*, *hlyJ*, *hlyK*, *hlyL*, *hlyM*, *hlyN*, *hlyO*, *hlyP*, *hlyQ*, *hlyR*, *hlyS*, *hlyT*, *hlyU*, *hlyV*, *hlyW*, *hlyX*, *hlyY*, *hlyZ*, *hlyA*, *hlyB*, *hlyC*, *hlyD*, *hlyE*, *hlyF*, *hlyG*, *hlyH*, *hlyI*, *hlyJ*, *hlyK*, *hlyL*, *hlyM*, *hlyN*, *hlyO*, *hlyP*, *hlyQ*, *hlyR*, *hlyS*, *hlyT*, *hlyU*, *hlyV*, *hlyW*, *hlyX*, *hlyY*, *hlyZ*, *hlyA*, *hlyB*, *hlyC*, *hlyD*, *hlyE*, *hlyF*, *hlyG*, *hlyH*, *hlyI*, *hlyJ*, *hlyK*, *hlyL*, *hlyM*, *hlyN*, *hlyO*, *hlyP*, *hlyQ*, *hlyR*, *hlyS*, *hlyT*, *hlyU*, *hlyV*, *hlyW*, *hlyX*, *hlyY*, *hlyZ*, *hlyA*, *hlyB*, *hlyC*, *hlyD*, *hlyE*, *hlyF*, *hlyG*, *hlyH*, *hlyI*, *hlyJ*, *hlyK*, *hlyL*, *hlyM*, *hlyN*, *hlyO*, *hlyP*, *hlyQ*, *hlyR*, *hlyS*, *hlyT*, *hlyU*, *hlyV*, *hlyW*, *hlyX*, *hlyY*, *hlyZ*, *hlyA*, *hlyB*, *hlyC*, *hlyD*, *hlyE*, *hlyF*, *hlyG*, *hlyH*, *hlyI*, *hlyJ*, *hlyK*, *hlyL*, *hlyM*, *hlyN*, *hlyO*, *hlyP*, *hlyQ*, *hlyR*, *hlyS*, *hlyT*, *hlyU*, *hlyV*, *hlyW*, *hlyX*, *hlyY*, *hlyZ*, *hlyA*, *hlyB*, *hlyC*, *hlyD*, *hlyE*, *hlyF*, *hlyG*, *hlyH*, *hlyI*, *hlyJ*, *hlyK*, *hlyL*, *hlyM*, *hlyN*, *hlyO*, *hlyP*, *hlyQ*, *hlyR*, *hlyS*, *hlyT*, *hlyU*, *hlyV*, *hlyW*, *hlyX*, *hlyY*, *hlyZ*, *hlyA*, *hlyB*, *hlyC*, *hlyD*, *hlyE*, *hlyF*, *hlyG*, *hlyH*, *hlyI*, *hlyJ*, *hlyK*, *hlyL*, *hlyM*, *hlyN*, *hlyO*, *hlyP*, *hlyQ*, *hlyR*, *hlyS*, *hlyT*, *hlyU*, *hlyV*, *hlyW*, *hlyX*, *hlyY*, *hlyZ*, *hlyA*, *hlyB*, *hlyC*, *hlyD*, *hlyE*, *hlyF*, *hlyG*, *hlyH*, *hlyI*, *hlyJ*, *hlyK*, *hlyL*, *hlyM*, *hlyN*, *hlyO*, *hlyP*, *hlyQ*, *hlyR*, *hlyS*, *hlyT*, *hlyU*, *hlyV*, *hlyW*, *hlyX*, *hlyY*, *hlyZ*, *hlyA*, *hlyB*, *hlyC*, *hlyD*, *hlyE*, *hlyF*, *hlyG*, *hlyH*, *hlyI*, *hlyJ*

SOLID™ next generation sequencing is an incredibly powerful tool that allows researchers to analyze a vast amount of genomic data in a relatively quick amount of time. Using this technology, we were able to sequence in one experimental run the whole genomes from 16 different *Salmonella* isolates representing 13 separate serogroups. With this data, it was possible to look at the evolutionary distance of multiple genomic markers, such as SPIs, across a broad spectrum of strains all at once. Previously, this type of in-depth analysis across a plethora of serogroups would not be possible in a timely manner. As part of the analyses, we were able to identify both large (such as loss of an ORF) and small (nucleotide sequence substitutions) genetic variations that may effect SPI function and possibly cause changes in disease severity and/or host specificity.

日期	星期	上午	下午	晚	备注
10月1日	星期一	语文	数学	英语	
10月2日	星期二	语文	数学	英语	
10月3日	星期三	语文	数学	英语	
10月4日	星期四	语文	数学	英语	
10月5日	星期五	语文	数学	英语	
10月6日	星期六	语文	数学	英语	
10月7日	星期日	语文	数学	英语	
10月8日	星期一	语文	数学	英语	
10月9日	星期二	语文	数学	英语	
10月10日	星期三	语文	数学	英语	
10月11日	星期四	语文	数学	英语	
10月12日	星期五	语文	数学	英语	
10月13日	星期六	语文	数学	英语	
10月14日	星期日	语文	数学	英语	
10月15日	星期一	语文	数学	英语	
10月16日	星期二	语文	数学	英语	
10月17日	星期三	语文	数学	英语	
10月18日	星期四	语文	数学	英语	
10月19日	星期五	语文	数学	英语	
10月20日	星期六	语文	数学	英语	
10月21日	星期日	语文	数学	英语	
10月22日	星期一	语文	数学	英语	
10月23日	星期二	语文	数学	英语	
10月24日	星期三	语文	数学	英语	
10月25日	星期四	语文	数学	英语	
10月26日	星期五	语文	数学	英语	
10月27日	星期六	语文	数学	英语	
10月28日	星期日	语文	数学	英语	
10月29日	星期一	语文	数学	英语	
10月30日	星期二	语文	数学	英语	
10月31日	星期三	语文	数学	英语	

Fig3A. Legend: Analysis of the protein identity for each CDS in SPI-1 was performed and color coded. Genes reported confirmed to be important for pathogenicity/virulence are highlighted in red in LT2 column. The average identity and average deviation from mean were calculated for each ORF across all isolates (minus isolates where CDS was reported not present). CDSs with the lowest % identity to reference sequence (red) and highest avg deviation (green) are shown. Isolates were ordered by the average identity of each ORF across the whole isolate. This isolate was also calculated for each isolate (minus ORFs 41-47) and listed at the bottom of each column. Isolates were positioned from left to right in order of highest to lowest % identity across the entire isolate.

Results: The analysis shows that the SPI-1 region in the Baldon isolate has diverged the least from the LT2 reference SPI-1 (99.62% identity), while the SPI-3 region in the Mississippi isolate has diverged the most (96.42% identity). The SPI-1 region in the Mississippi isolate has diverged the least (99.62% identity) from the *o/crc* and *spsB*. Interestingly, both *ygbA* and *orcF* have been shown NOT to be important for virulence in LT2, while *avrA*, CDSs with the highest protein sequence variation include *ygbA*, *prgI* and *spsD*. CDSs corresponding to *STM2901* through *STM2908* are all missing in at least 1 isolate, and may suggest that they were inherited in the same horizontal acquisition event as the rest of the island, but are not important for pathogenicity and so are not listed.

Island	SP1	SP2	SP3	SP4	SP5	SP6	SP7	SP8	SP9	SP10	SP11	SP12	SP13	SP14 (avg SP1-2)		
Genes missing in at least 1 isolate	<i>avrA</i>	<i>ssaA</i>	STM752	STM4205	STM1089	STM2744	STM2952	all	all	STY2878	all	SC1240	SC1249	ita	STM317	all
	STM2901	<i>supR</i>			STM1093	STM2075	STM2903					<i>scgB</i>	SC1250	vrf	STM318	
	STM2902		STM754			STM2276	STM2904					SC1242	capH	CG2248	STM319	
	STM2903		<i>muhI</i>			STM2825	STM2904.IN					<i>mnpA</i>	SC1240		STM320	
	STM2904					STM3287	STM2905					page			STM321	
	STM2905					STM2386	STM2907						SC1245		STM322	
	<i>gphH</i>					STM2909							SC1246		STM323	
	STM2908					STM2901							<i>oriA</i>			

Fig 3B. Legend: The analysis shown in figure 3A was performed on the multiple SPIs. Orthologues not present in at least 1 isolate in this study are listed here. Genes known to be required for pathogenesis in the reference strain are highlighted in red.

[illegible]

Fig4. Legend: Phylogenetic analysis was performed on 10 SPI orthologues scattered throughout SPIs 1, 2 and 5. Pairwise distance calculations (averages listed in fig4b) and phylogeny reconstruction using a neighbouring-joining algorithm (averages listed in fig4c) for *saeC* shown in fig4d) were performed. Cluster Analysis of the phylogeny reconstruction was used to determine which isolates cluster together. Clusters were numbered according to distance from reference isolate in the tree (see table 1).
Results: From the analysis, we were able to make 3 observations. 1. Similar gene cluster patterns can be found for several of the CDs and suggests that many of these CDs have co-evolved (such as *saeCDP* and *saeC*) in isolates. Also, the *saeC* gene is conserved in all isolates. 2. The *saeC* gene is conserved in all Senftenberg isolates have diverged in their own cluster from the other *saeC* genes. Since both of these strains are derived from a bovine source, it may suggest that alterations in *Sae* (helping to determine host specificity, 3. Isolates Uganda and Kenya have evolved independently, as demonstrated by an almost total lack of common clusters (only shared clusters suggesting part of the same group) (E).

- SPIs 1-5, 9, and 13 are universal among a wide range of serogroups.
- SPIs 10, 15, 17, and 19-21 seem to have a highly serovar-restricted profile.
- SPIs 7 and 11 seemed conserved in multiple serogroups but located in different genomic regions.
- Most Prophage insertions are random and are not conserved between strains.
- ORF analysis for each SPI may be helpful in re-defining pathogenicity island boundaries and in determining if a virulence factor has a serovar-restricted or a more universal phenotype
- Phylogenetic analysis of SPI CDSs may help determine which genes may be shaping disease outcome and host specificity by looking for individual or groups of virulence factors that have evolved together.