

GeneBridge 4 Human/Hamster Radiation Hybrid Panel

Description

The GeneBridge 4 Human/Hamster Radiation Hybrid Panel consists of low resolution 93 Radiation Hybrid clones of the whole human genome. This is a subset of the 199 clone panel developed in collaboration between the labs of Peter Goodfellow and Jean Weissenbach (Walter, M. A., Spillett, D. J., Thomas, P., Weissenbach, J., and Goodfellow, P. N. (1994). A method for constructing radiation hybrid maps of whole genomes. *Nature Genetics* 7: 22 - 28). A human cell line (donor, high frequency lysogeny (HFL)) was exposed to 3,000 rad of X-rays and then fused with thymidine-deficient hamster recipient cells (A23), creating a panel of hybrids with 1,000 kb resolution.

Contents

The GeneBridge 4 Radiation Hybrid Panel contains a vial of DNA from each of 93 radiation hybrid clones, plus the two control DNAs (HFL donor and A23 recipient, 5 µg each). Each vial contains 25 ng / µl of DNA in Tris-EDTA buffer.

Perform PCR using 25 ng of DNA in 10 µl PCR reactions.

Catalog nos.	Amount
RH02.02	2 µg of each hybrid
RH02.05	5 µg of each hybrid
RH02.50	50 µg of each hybrid

Server for Mapping STSs

<http://www-genome.wi.mit.edu/> is a publicly available server that allows mapping sequence tagged sites (STSs) relative to the Whitehead Institute / MIT Center for Genome Research's radiation hybrid map of the human genome, which is constructed with the GeneBridge 4 Radiation Hybrid Panel.

Hybrid Order for Entry in the Server

On the next page, these hybrids are listed in the order in which they must be entered in the server. (For reading convenience only, we group hybrids in sets of five.)

Storage

Store the Radiation Hybrid DNA preparations at -20 °C.

©2001-2002 Invitrogen Corporation. All rights reserved.

25-0549B 062502

This product is distributed for laboratory research use only. CAUTION: Not for diagnostic use. The safety and efficacy of this product in diagnostic or other clinical uses has not been established.

For technical questions about this product, call the Invitrogen TECH-LINESM 800 955 6288

RH PCR and Scoring - GeneBridge 4 Panel

Project:

Sequence tagged site (STS) name:

Temperature / buffer:

Fragment length:

Chromosome:

Date of primary PCR run:

Date of gel running:

Scored by:

HFL (human)

A23 (hamster)

A4 - A5 - AA5 - AA7 - B2

B3 - B9 - BB1 - BB6 - BB8

BB10 - BB12 - C3 - C11 - CC8

D1 - D7 - DD2 - DD5 - DD8

E2 - E4 - E6 - E11 - F6

F7 - F13 - G1 - G5 - G6

G7 - G11 - H1 - H8 - H9

H12 - I1 - I4 - J2 - J5

J9 - K5 - K7 - K8 - K9

K12 - L3 - L4 - L6 - M4

M5 - N3 - N5 - N6 - N7

N12 - O5 - O10 - P2 - P9

P11 - Q2 - Q4 - R1 - R2

R3 - R5 - R6 - R10 - R12

S3 - S6 - S10 - S12 - T3

T4 - T10 - T11 - U1 - U3

V2 - V3 - V7 - V8 - W1

Y4 - Y8 - Y9 - Z5 - Z6

Z9 - Z11 - Z12