

DS-32 Matrix Standard Kit (Dye Set F)

SeqStudio™ Flex, SeqStudio™, 3500, and 3130 series instruments

Catalog Number 4345831

Pub. No. 4362855 Rev. E

 **WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from thermofisher.com/support.

Product description

The DS-32 Matrix Standard Kit (Dye Set F) is used to perform spectral calibrations when analyzing DNA fragments labeled with 5-FAM™, JOE™, NED™, and ROX™ dyes. (The ROX™ dye is used to label the size standard.) The matrix standard contains four DNA fragments. Each fragment is labeled with a different dye from the dye set.

For more information on spectral calibration, see the *DNA Fragment Analysis by Capillary Electrophoresis User Guide* (Pub. No. 4474504).

Contents and storage

Contents	Amount	Storage
DS-32 Matrix Standard in TE buffer	1 tube	Store at 2–8°C, protected from light. ^[1] Do not freeze.

^[1] The kit is stable for 1 year when stored at 2–8°C.

Required materials not supplied

Item	Cat. No.	
Hi-Di™ Formamide	4311320	
MicroAmp™ Fast Optical 96-Well Reaction Plate, 0.1 mL	4346907	
MicroAmp™ Optical 96-Well Reaction Plate	N8010560	
MicroAmp™ Optical 384-Well Reaction Plate	4343370	
Septa		
SeqStudio™ Flex and 3500 series	8-Strip Septa 3500/Flex Series (Qty 24) 96-Well Septa 3500/Flex Series	4410701 4412614
SeqStudio™	384-Well Septa 3500/Flex Series	4412520
SeqStudio™	Septa for SeqStudio™ Genetic Analyzer, 96 well	A35641
3130 series	Plate Septa, 96 well	4315933

Guidelines for use

- For more information on the use of matrix standards, see the instrument user guide or getting started guide.
- To prepare the matrix standard dilution, combine the appropriate volumes of matrix standard and Hi-Di™ Formamide (Cat. No. 4311320). Dilution volumes vary depending on the instrument.
- Use the matrix standard within 2 hours of preparation.
- Do not add size standard to the matrix standard.
- Discard any unused reagent that has been diluted in Hi-Di™ Formamide.

Prepare the standard

1. Vortex the matrix standard tube for 5–10 seconds to mix, then centrifuge for 3–5 seconds to bring the mixture to the bottom and eliminate air bubbles.
2. Combine the volumes of matrix standard and Hi-Di™ Formamide appropriate for the instrument. See “Component volumes and well location for the prepared standard” on page 2.
3. Vortex for 5–10 seconds, then centrifuge for 3–5 seconds.
4. Dispense the prepared standard into the appropriate wells of a reaction plate. See “Component volumes and well location for the prepared standard” on page 2.
5. Cover the plate with adhesive film, then centrifuge for 3–5 seconds.
6. Denature the DNA fragments:
 - a. Incubate the mixture at 95°C for 5 minutes.
 - b. Incubate the mixture at 4°C, or on ice, for ≥2 minutes.
7. Remove the adhesive film, then cover the plate with septa.
8. Centrifuge for 3–5 seconds.
9. Assemble the plate with the retainer and base, then load on the instrument.

Note: The SeqStudio™ Genetic Analyzer does not require a retainer and base.
10. Immediately perform the spectral calibration.

For information on setting up the run, see the instrument user guide.

Component volumes and well location for the prepared standard

Table 1 SeqStudio™ Flex Series Genetic Analyzer

Component	Volume		Well location for the prepared standard	
	8-capillary array	24-capillary array	96-well plate	384-well plate
DS-32 Matrix Standard	2 µL	6 µL	Dispense 10 µL of the prepared standard into wells of a 96-well plate: <ul style="list-style-type: none"> • 8-capillary array—8 wells (for example, A1–H1) • 24-capillary array—24 wells (for example, A1–H3, A4–H6, A7–H9, or A10–H12) 	Dispense 5 µL of the prepared standard into wells of a 384-well plate: <ul style="list-style-type: none"> • 24-capillary array—24 wells (for example, A1, A3, A5; C1, C3, C5; E1, E3, E5; G1, G3, G5; I1, I3, I5; K1, K3, K5; M1, M3, M5; O1, O3, O5)
Hi-Di™ Formamide	123 µL	369 µL		
Total volume	125 µL	375 µL		

Table 2 SeqStudio™ Genetic Analyzer

Component	Volume	Well location for the prepared standard
	4-capillary array	
DS-32 Matrix Standard	1 µL	Dispense 10 µL of the prepared standard into wells of a 96-well plate: <ul style="list-style-type: none"> • 4 wells (for example, A1–D1)
Hi-Di™ Formamide	49 µL	
Total volume	50 µL	

Table 3 3500/3500xL Genetic Analyzer

Component	Volume		Well location for the prepared standard
	8-capillary array	24-capillary array	
DS-32 Matrix Standard	4 µL	4 µL	Data Collection Software v3 and later: Dispense 10 µL of the prepared standard into wells of a 96-well plate: <ul style="list-style-type: none"> • 8-capillary array—8 wells (for example, A1–H1) • 24-capillary array—24 wells (for example, A1–H3, A4–H6, A7–H9, or A10–H12) Note: If you place the standard in wells that do not correspond to injection position 1, specify the starting well position in the software. Data Collection Software v1, v1.1, and v2: Dispense 10 µL of the prepared standard into wells of a 96-well plate: <ul style="list-style-type: none"> • 8-capillary array—8 wells: A1–H1 • 24-capillary array—24 wells: A1–H3
Hi-Di™ Formamide	246 µL	246 µL	
Total volume	250 µL	250 µL	

Table 4 3130/3130x/ Genetic Analyzer

Component	Volume		Well location for the prepared standard
	36-cm array	50-cm array	
DS-32 Matrix Standard	10 µL	5 µL	Dispense 10 µL of the prepared standard into wells of a 96-well plate: <ul style="list-style-type: none"> • 16-capillary array—16 wells: A1–H2 • 4-capillary array—4 wells: A1–D1
Hi-Di™ Formamide	190 µL	195 µL	
Total volume	200 µL	200 µL	

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Thermo Fisher Scientific Baltics UAB | V.A. Graiciuno 8, LT-02241 | Vilnius, Lithuania
 For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision history: Pub. No. 4362855

Revision	Date	Description
E	31 January 2022	Added information for the SeqStudio™ Flex Series Genetic Analyzer. Added required materials table.
D	10 April 2020	Added new formulation for the SeqStudio™ Genetic Analyzer. Removed 3100 series instruments. Added TE buffer to the contents description. Added vortex and centrifuge times. Added information for Data Collection Software v1, v1.1, and v2. Changed the manufacturing address to Vilnius. Updated format and licensing.
C	30 June 2015	Baseline for this revision.

Important Licensing Information: This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2022 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

