

Agilent Seahorse XF HS Mini Analyzer Consumables



For use with Seahorse XF HS Mini Analyzer model number series: S7852A

FluxPaks

Part No.	Description	No. Cartridges/Miniplates
103723-100	XF HS Mini FluxPak	12/12
103724-100	XF HS Mini FluxPak (PDL plates)	12/12
103721-100	XFp FluxPak (PDL plates)	12/12
103022-100	XFp FluxPak	12/12

Miniplates and Miniplate Carriers

Part No.	Description	No. Miniplates or Trays
103725-100	XF HS Miniplates	12
103727-100	XF HS PDL Miniplates	12
103722-100	XFp PDL Miniplates	12
103025-100	XFp Cell Culture Miniplates	12
103057-100	XFp Carrier Tray	4

Assay Kits and Reagents

Part No.	Description	No. of 8 Well Assays
103591-100	Seahorse XFp Real-Time ATP Rate Assay Kit	6
103010-100	Seahorse XFp Cell Mito Stress Test Kit	6
103346-100	Seahorse XFp Glycolytic Rate Assay Kit	6
103017-100	Seahorse XFp Glycolysis Stress Test Kit	6
103275-100	Seahorse XFp Cell Energy Phenotype Test Kit	12
103270-100	Seahorse XFp Mito Fuel Flex Test Kit	6
102504-100	Seahorse XF Plasma Membrane Permeabilizer	60
102720-100	Seahorse XF Palmitate-BSA FAO Substrate	30

Starter Packs

Part No.	Description	No. of 8 Well Assays
103707-100	Seahorse XFp Cell Mito Stress Test Starter Pack: Includes one XFp Cell Mito Stress Test Kit (103010-100), 500 mL XF DMEM medium pH 7.4 (103575-100), 50 mL XF 1.0 M Glucose Solution (103577-100), 50 mL XF 100 mM Pyruvate Solution (103578-100), and XF 200 mM Glutamine Solution (103579-100).	6
103709-100	Seahorse XF Glycolytic Rate Assay Starter Pack: Includes one XFp Glycolysis Rate Assay kit (103346-100), 500 mL XF DMEM medium pH 7.4 (103575-100), 50 mL XF 1.0 M Glucose Solution (103577-100), 50 mL XF 100 mM Pyruvate Solution (103578-100), and XF 200 mM Glutamine Solution (103579-100).	6
103676-100	Seahorse XFp Real-Time ATP Rate Assay Starter Pack: Includes one XFp Real-Time ATP Rate Assay Kit (103591-100), 500 mL XF DMEM medium pH 7.4 (103575-100), 50 mL XF 1.0 M Glucose Solution (103577-100), 50 mL XF 100 mM Pyruvate Solution (103578-100) and XF 200 mM Glutamine Solution (103579-100).	6

Assay Media and More

Part No.	Description	Quantity
103680-100	Seahorse XF DMEM assay medium pack: Includes 500 mL XF DMEM medium pH 7.4 (103575-100), 50 mL XF 1.0 M Glucose Solution (103577-100), 50 mL XF 100 mM Pyruvate Solution (103578-100), and XF 200 mM Glutamine Solution (103579-100).	
103681-100	Seahorse XF RPMI assay medium pack: Includes 500 mL XF RPMI medium pH 7.4 (103576-100), 50 mL XF 1.0 M Glucose Solution (103577-100), 50 mL XF 100 mM Pyruvate Solution (103578-100), and XF 200 mM Glutamine Solution (103579-100).	
103575-100	Seahorse XF DMEM medium, pH 7.4	500 mL
103576-100	Seahorse XF RPMI medium, pH 7.4	500 mL
103577-100	Seahorse XF 1.0 M Glucose Solution	50 mL
103578-100	Seahorse XF 100 mM Pyruvate Solution	50 mL
103579-100	Seahorse XF 200 mM Glutamine Solution	50 mL
103193-100	Seahorse XF Base Medium	100 mL
103334-100	Seahorse XF Base Medium	500 mL
102353-100	Seahorse XF Base Medium	2 x 1L
103335-100	Seahorse XF Base Medium, without phenol red	500 mL
100840-000	Seahorse XF Calibrant Solution	500 mL
103059-000	Seahorse XF Calibrant Solution	100 mL

www.agilent.com

For Research Use Only. Not for use in diagnostic procedures.

This information is subject to change without notice.

© Agilent Technologies, Inc. 2019, 2020
Printed in the USA, January 24, 2020
5991-8589EN