

Maxwell® 16 Polyhistidine Protein Purification Kit

Learn more. Meet Maxwell® 16.

The Maxwell® 16 Polyhistidine Protein Purification Kit is used with the Maxwell® 16 Instrument to provide an easy method for the efficient, automated purification of polyhistidine-tagged protein from bacterial cultures and other sample types including mammalian and insect cells. With minor modifications, the reagents can also be used for purification of HQ-tagged proteins from bacterial cultures.

The Maxwell® Instrument is supplied with a preprogrammed purification procedure and reagent cartridges specifically designed to maximize simplicity and con...

Expand to Read More »



Prices valid for US customers only

PRODUCT	SIZE	CONC.	CATALOG #	*LIST PRICE	ORDER QTY	
Maxwell® 16 Polyhistidine Protein Purification Kit	48 preps	-	AS1060	\$ 332.00		Add to cart
Protocol MSDS Components						

Available Separately

PRODUCT	SIZE	CONC	CATALOG #	*LIST PRICE	ORDER QTY	.
SEV Plungers Components	50/pk	-	AS5201	\$ 20.00		Add to cart
SEV Elution Tubes Components	50/pk	-	AS5101	\$ 6.00		Add to cart

You Also May Be Interested In



Maxwell® 16 Blood DNA Purification Kit

Maxwell® 16 Cell DNA Purification Kit

Maxwell® 16 Tissue DNA Purification Kit

Maxwell® 16 Total RNA Purification Kit

Storage Conditions

Store at 4°C.

For product intended use please see Patents & Disclaimers tab.

Maxwell® 16 Intra-Run and Inter-Run Consistency

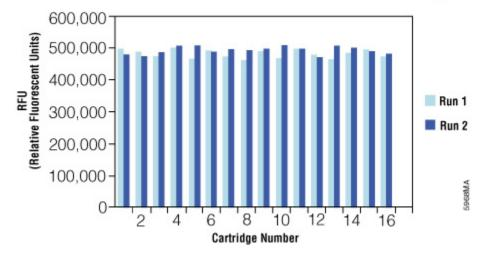


Figure 1. Intra-run and inter-run consistency of luciferase purification using the Maxwell® 16 Polyhistidine Protein Purification Kit and the Maxwell® 16 Instrument.

Protocols

Maxwell® 16 Polyhistidine Protein Purification Kit Technical Manual

Articles

Maxwell® 16 Polyhistidine Protein Purification Kit

Related Products

DNA IQ™ Reference Sample Kit for Maxwell® 16

Maxwell® 16 System DNA Purification Kits

Maxwell® 16 System RNA Purification Kits

Maxwell® 16 Instrument for Nucleic Acid and Protein Purification

Related Citations

Small-scale, semi-automated purification of eukaryotic proteins for structure determination. 2008 *J. Struct. Funct. Genomics* **8**, 153-166.

Service and Support

Service and Support

Maxwell® 16 System Firmware

Use Restrictions

AS1060, AS5201, AS5101 For Research Use Only. Not for Use in Diagnostic Procedures.

Patents - Disclaimers

AS1060 U.S. Pat. Nos. 7,112,552 and 7,354,750.

AS1060 This product is licensed for use under U.S. Pat. No. 6,174,704.

View All Reviews »

Review This Product »

by