



# 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	QTY	
Maxwell® 16 Polyhistidine Protein Purification Kit	48 preps	-	AS1060	\$ 345.00	<input type="text"/>	<a href="#">Add to cart</a>
<a href="#">Protocol</a>						
<a href="#">MSDS</a>						
<a href="#">Components</a>						

## Available Separately

PRODUCT	SIZE	CONC.	CATALOG #	*LIST PRICE	QTY	
SEV Plungers	50/pk	-	AS5201	\$ 35.00	<input type="text"/>	<a href="#">Add to cart</a>
<a href="#">Components</a>						
SEV Elution Tubes	50/pk	-	AS5101	\$ 36.00	<input type="text"/>	<a href="#">Add to cart</a>
<a href="#">Components</a>						

## Storage Conditions

Store at 4°C.

For product intended use please see Patents & Disclaimers tab.

## 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**