

## MsrB2 protein (Human)

aa: 42-182, 6xHis

Store at 4°C (desiccated)

**Catalog #** MB201  
MB201-XL

**Product Description:** Human MSRB2 protein (MB201) [accession # Q9Y3D2 (MSRB2\_HUMAN)] has been produced and purified from a bacterial expression system. The recombinant protein is N-terminally 6x Histidine tagged and is comprised of amino acids 42-182 [5]. MB201 has an approximate molecular weight of 17 kDa.

**Amount:** MB201 50 µg  
MB201-XL 1 mg

**Lot:** 015

TEST	SPECIFICATION	LOT RESULTS
<b>Appearance</b>	White lyophilized powder	White lyophilized powder
<b>Purity by Coomassie Gel</b>	>90%	>90%
<b>Protein Quantitation with Precision Red Protein Assay</b>	MB201 50-60µg	53 µg
<b>Reagent (Cytoskeleton Cat. # ADV02)</b>	MB201-XL 1.0-1.2 mg	1.13 mg
<b>Biological Activity Assay #1: Subtilisin cleavage</b>	<10% cleavage of Mical1 oxidized actin. >90% cleavage of Mical1 oxidized actin treated with MsrB2	<10% >90%
<b>Biological Activity Assay #2: Pyrene actin polymerization assay</b>	MsrB2 (1.2 µM) treatment of oxidized pyrene actin (0.05 mg/ml) is able to restore polymerization activity from 0-5% to 80-90% the RFU values of native pyrene actin.	Oxidized actin 0% MsrB2 treated actin 90%
<b>QC Release Date</b>	May 8 <sup>th</sup> 2019	



Kim Middleton  
QC Manager