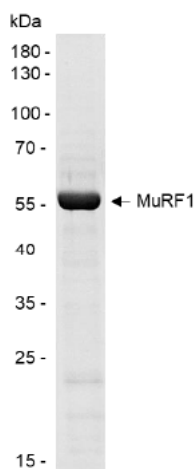


## Product Information

<b>Description:</b>	Human MuRF1 recombinant protein (Genbank Accession No. NP_115977), full length with N-terminal SUMO- and His-tags, expressed in <i>E. coli</i> .
<b>Background:</b>	MuRF1 (Muscle-specific RING-finger protein 1) is an E3 (ubiquitin protein ligase) that is involved in the conjugation of ubiquitin to target substrates along with E1 and E2 enzymes. MuRF1 is also known as TRIM63 (tripartite motif containing 63) containing a RING-finger/B-box/coiled-coil tripartite fold and is a member of the RING-domain E3 family. MuRF1 has been implicated as a component of atrophy-associated accelerated proteolysis. It is upregulated in a number of different models of atrophy and interacts with numerous myofibrillar proteins.
<b>Species:</b>	Human
<b>Tag:</b>	His-SUMO
<b>Concentration:</b>	0.40 mg/ml
<b>Expression System:</b>	<i>E. coli</i>
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	50 mM Tris Buffer, pH 7.5, 150 mM NaCl, 10% glycerol, 5 mM DTT
<b>MW:</b>	55 kDa
<b>Genbank Accession:</b>	NP_115977
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
<b>Biological Activity</b>	Typical enzyme concentration of 100 nM - 5 mM is used for in vitro conjugation, depending on conditions.
<b>Assay Conditions:</b>	3.125 nM MuRF1 was tested in a TR-FRET assay showing robust E3 ligase activity.
<b>Applications:</b>	Useful as an E3 ligase in ubiquitin-conjugating reactions and to analyze upregulation/downregulation of E3 enzymes.

## Quality Control Data

## 10-20% SDS-PAGE Coomassie Staining



## MuRF1 Activity

