

# MuRF1 (Muscle-specific RING finger protein 1)

Cat. # UB301

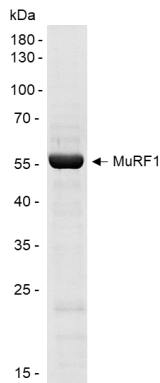
## Background

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 several different models of atrophy and interacts with numerous myofibrillar proteins.

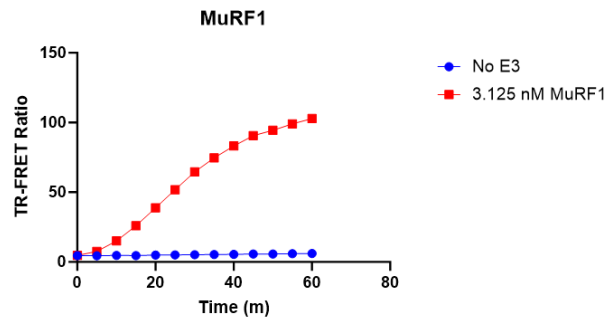
## Product Information

<b>Purity</b>	≥ 90% by SDS-PAGE
<b>Molecular Weight</b>	40 kDa (Without Tag), ~55 kDa with SUMO tag
<b>Quantity</b>	25 µg
<b>Physical State</b>	Liquid
<b>Species</b>	Human
<b>Source</b>	<i>E. coli</i>
<b>Tag</b>	His6-SUMO
<b>Activity</b>	Typical enzyme concentration of 100 nM - 5 mM is used for in vitro conjugation, depending on conditions.
<b>Storage</b>	-80° C. Avoid repeated freeze/thaw cycles

## Product QC



**SDS-Page Analysis of purified MuRF1.** Two µg of the protein was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



**Activity Assay of MuRF1.** 3.125 nM MuRF1 was tested in a TR-FRET assay showing robust E3 ligase activity.

## References

- Adams, V., et al., J Cachexia Sarcopenia Muscle., 2022. 13(3):1565-1581.
- Baehr, LM., et al., Function (Oxf)., 2021. 2(4)zqab029.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic use.  
Copyright © 2009 LifeSensors, Inc. All Rights Reserved