

MuRF2 (Muscle-specific RING finger protein 2)

Cat. # UB305

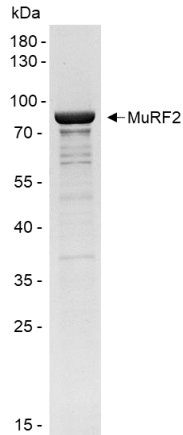
Background

MuRF2 is a RING domain E3 ligase that is involved in the conjugation of ubiquitin to target substrates. MuRF2 has been demonstrated to function with the E2 enzyme UBE2D3 (UbcH5c) in vitro. MuRF2 is also known as TRIM55 (tripartite motif-containing 55) containing a RING-finger/B-box/coiled-coil tripartite fold. MuRF2 has been implicated along with MuRF1 as regulators of protein degradation in striated muscle.

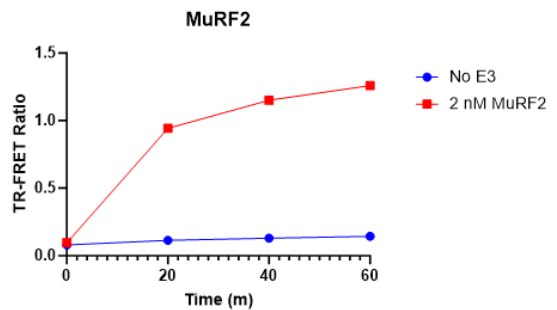
Product Information

Purity	≥ 85% by SDS-PAGE
Molecular Weight	72.7 kDa (with tag), 60.5 kDa (without tag)
Quantity	25 µg
Physical State	Liquid
Species	Human
Source	<i>E. coli</i>
Tag	His6-SUMO
Activity	Typical enzyme concentration of 100 nM - 5 mM is used for in vitro conjugation, depending on conditions.
Storage	-80° C. Avoid repeated freeze/thaw cycles

Product QC



SDS-Page Analysis of purified MuRF2. Two µg of the protein was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



Activity Assay of MuRF2. 2 nM MuRF2 was tested in a TR-FRET assay showing robust E3 ligase activity.

References

1. Bian, H., et al., FEBS Open Bio., 2018. 8(2):234-243.
2. Silvestre, JG., et al., Braz J Med Biol Res., 2019. 52(9):e8551.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic use.

Copyright © 2009 LifeSensors, Inc. All Rights Reserved