

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/Bbox/coiled-coil tripartite fold. MuRF2 has been implicated along with MuRF1 as regulators of protein degradation in striated muscle.

Alternate Names

Tripartite Motif-Containing Protein 55 (TRIM55), Muscle-Specific RING Finger Protein 2, Ring Finger Protein 29 (RNF29)

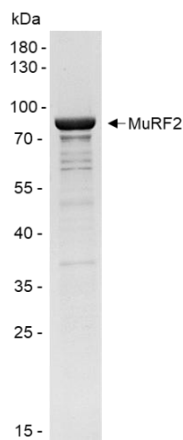
Application(s)

In vitro conjugation assay

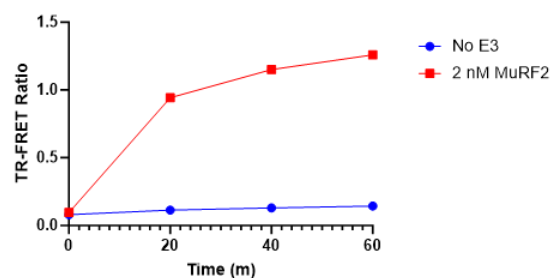
Product Specifications

Affinity tag	His6-SUMO
Purity	> 85% by SDS-PAGE
Molecular Weight	72.7 kDa (with tag), 60.5 kDa (without tag)
Quantity	25 µg
Species	Human
Expression System	E. coli
Physical State	Liquid
Buffer	50 mM Tris, 150 mM NaCl, 5 mM DTT (fresh), 10% glycerol
Activity	A typical enzyme concentration of 1-100 nM is used for in vitro conjugation, depending on experimental conditions.
Storage	Store at -80°C. Avoid repeated freeze/thaw cycles

Product QC



MuRF2



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 uses
Copyright © 2025 LifeSensors, Inc. All Rights Reserved