

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.

Alternate names: E3 ubiquitin-protein ligase TRIM55, MURF2, MuRF-2, MURF-2, Muscle-specific RING finger protein 2, RING finger protein 29, RNF29, Tripartite motif-containing protein 55.

Product Information

Purity:	≥ 95% by SDS-PAGE
Molecular Weight:	72.7kDa (with tag)
Quantity:	25 µg
Physical State:	Liquid
Source:	Recombinant Human
Tag:	His ₆ SUMO
Activity:	Typical enzyme concentration 20nM -1 mM is used for <i>in vitro</i> conjugation depending on assay conditions
Storage:	-80° C. Avoid repeated freeze/thaw cycles

References

1. Willis, M.S., C. Ike, L. Li, D.Z. Wang, D.J. Glass, and C. Patterson, Muscle ring finger 1, but not muscle ring finger 2, regulates cardiac hypertrophy *in vivo*. *Circ Res*, 2007. **100**:456-9.
2. Witt, S.H., H. Granzier, C.C. Witt, and S. Labeit, MURF-1 and MURF-2 target a specific subset of myofibrillar proteins redundantly: towards understanding MURF-dependent muscle ubiquitination. *J Mol Biol*, 2005. **350**:713-722.
3. Witt, C.C., S.H. Witt, S. Lerche, D. Labeit, W. Back, and S. Labeit, Cooperative control of striated muscle mass and metabolism by MuRF1 and MuRF2. *EMBO J*, 2008. **27**:350-360.
4. Moriscot, A.S., I.L. Baptista, J. Bogomolovas, C. Krohne, S. Hirner, H. Granzier, and S. Labeit, MuRF1 is a muscle fiber-type II associated factor and together with MuRF2 regulates type-II fiber trophicity and maintenance. *J Struct Biol*. 2010, **170**:344-53.

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