

MuRF3 (Muscle-specific RING finger protein 3)

Cat. # UB306

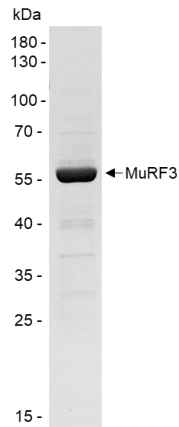
Background

MuRF3 is a RING domain E3 ligase that is involved in the conjugation of ubiquitin to target substrates. MuRF3 has been demonstrated to function with the E2 enzyme UBE2D3 (UbcH5c) in vitro. MuRF3 is also known as TRIM54 (tripartite motif-containing 54) containing a RING-finger/B-box/coiled-coil tripartite fold. MuRF3 has been implicated along with MuRF1 as a regulator of protein degradation in striated muscle. MuRF3 also plays an important role in maintaining cardiac function after myocardial infarction.

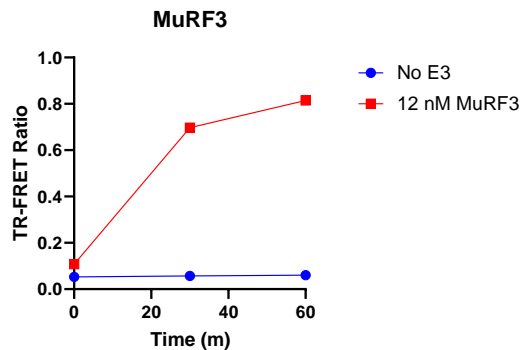
Product Information

Purity	≥ 90% by SDS-PAGE
Molecular Weight	52.4 kDa (with tag), 40.3 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 MuRF3. Two µg of the protein was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



Activity Assay of MuRF3. 12 nM MuRF3 was tested in a TR-FRET assay showing robust E3 ligase activity.

References

1. Cao, H., et al., Asia Pac J Clin Oncol., 2022. 18(6):669-677.
2. Zhu, J., et al., Front Oncol., 2021. 11:759842.

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

Copyright © 2009 LifeSensors, Inc. All Rights Reserved