

Background

The E3 ubiquitin ligase TRAF6, together with UBE2N and UBE2V1, mediates the formation of Lys63-linked polyubiquitin chains on proteins such as IKBKG, IRAK1, AKT1, and AKT2. TRAF6 also facilitates the ubiquitination of free (unanchored) polyubiquitin chains, which leads to the activation of MAP3K7. It plays a critical role in the development of functional osteoclasts. Additionally, TRAF6, along with MAP3K8, mediates CD40 signaling, which activates ERK in B cells and macrophages and may influence the regulation of immunoglobulin production.

Alternate Names

TNF Receptor Associated Factor 6, RING Finger Protein 85, RING-Type E3 Ubiquitin Transferase TRAF6, E3 Ubiquitin-Protein Ligase TRAF6, Interleukin-1 Signal Transducer

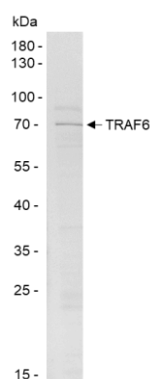
Application(s)

Ubiquitin ligation reactions

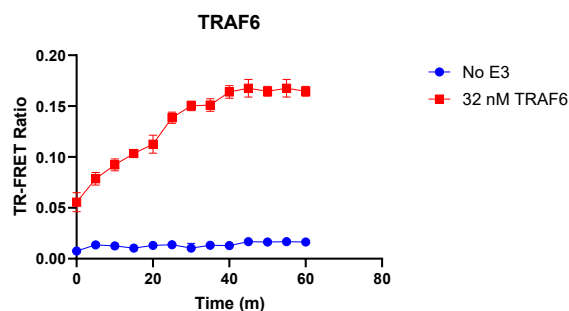
Product Specifications

Affinity tag	His ₆ -SUMO
Purity	> 85% by SDS-PAGE
Molecular Weight	70 kDa
Quantity	25 µg
Species	Human
Expression System	<i>E. coli</i>
Physical State	Liquid
Buffer	1X PBS, 10% glycerol
Activity	A typical enzyme concentration of 10-100 nM is used for in vitro conjugation, depending on assay conditions
Storage	Store at -80°C. Avoid repeated freeze/thaw cycles

Product QC



SDS-Page analysis of purified Traf6. Two µg of the protein was loaded onto a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



Activity Assay of Traf6. TRAF6 at a concentration of 32 nM was tested in a TR-FRET assay, demonstrating robust E3 ligase activity.

References

1. Min, Y., et al., Autophagy, 2018. 14(7):1347-1358.
2. Wang, YT., et al., Cell Rep, 2022 38(8):110354

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2025 LifeSensors, Inc. All Rights Reserved