

Traf6

Cat. # UB312

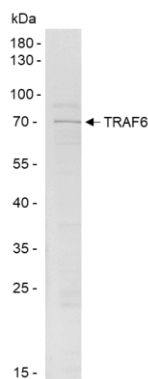
Background

TRAF6 is an E3 ubiquitin ligase that, together with UBE2N and UBE2V1, mediates the synthesis of 'Lys-63'-linked-polyubiquitin chains conjugated to proteins, such as IKK β , IRAK1, AKT1 and AKT2. Traf6 also mediates ubiquitination of free/unanchored polyubiquitin chain that leads to MAP3K7 activation. Traf6 may be essential for the formation of functional osteoclasts. Together with MAP3K8, Traf6 mediates CD40 signals that activate ERK in B-cells and macrophages, and thus may play a role in the regulation of immunoglobulin production.

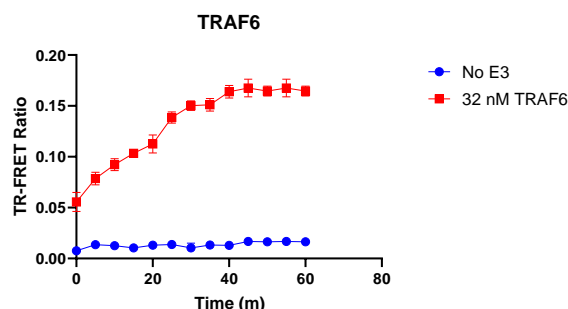
Product Information

Purity	≥ 85% by SDS-PAGE
Molecular Weight	70 kDa
Quantity	25 µg
Physical State	Liquid
Species	Human
Source	<i>E. coli</i>
Tag	His6-SUMO
Activity	Typical enzyme concentration 50nM -1 µM is used for in vitro conjugation depending on assay conditions
Storage	-80° C. Avoid repeated freeze/thaw cycles

Product QC



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



Activity Assay of Traf6. 32 nM Traf6 was tested in a TR-FRET assay showing a 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 use.
Copyright © 2013 LifeSensors, Inc. All Rights Reserved