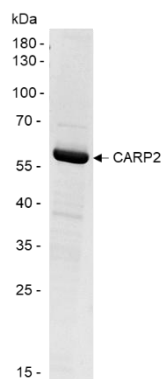


Background	CARP2 (caspase 8/10-associated RING protein 2) is a RING-domain E3 (ubiquitin protein ligase) involved in the conjugation of ubiquitin to target substrates, along with E1 and E2 enzymes. CARPs (CARP1 and CARP2, 77% identity) also belong to the IAP family (inhibitors of apoptosis proteins) inhibiting the activation of DED (death effector domain)-containing caspase proteins (8/10). In addition to targeting caspases 8/10, CARPs have been shown to target phosphorylated p53 for degradation in an Hdm2-independent manner. Furthermore, CARP2 has been shown to be a negative regulator of TNF-induced NF- κ B activation by targeting RIP for degradation.
Alternate Names	RING Finger And FYVE-Like Domain-Containing Protein 1, RING-Type E3 Ubiquitin Transferase Rififylin, RING Finger Protein 189 (RNF189), Caspase Regulator CARP2, Ring Finger And FYVE-Like Domain Containing 1, FYVE-RING Finger Protein SAKURA, RING Finger Protein 34-Like (RNF34L), Fring, Rififylin
Application(s)	In vitro conjugation assay

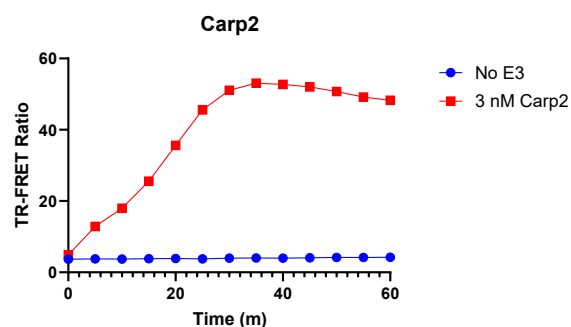
Product Specifications

Affinity tag	His6-SUMO
Purity	> 90% by SDS-PAGE
Molecular Weight	41 kDa (Without Tag), ~55 kDa with SUMO tag
Quantity	25 μ g
Species	Human
Expression System	<i>E. coli</i>
Physical State	Liquid
Buffer	50 mM Tris, 150 mM NaCl, 10 mM DTT
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



SDS-Page Analysis of purified CARP2. Two μ g of the protein was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



Activity Assay of Carp2. 3 nM Carp2 was tested in a TR-FRET assay showing robust E3 ligase activity.

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

References

1. Sharma, R., et al., FEBS J., 2023. 290(14):3580-3594.
2. Weng, X., et al., Int J Biol Macromol., 2023. 224:713-724.

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