

ITCH

Cat. # UB317

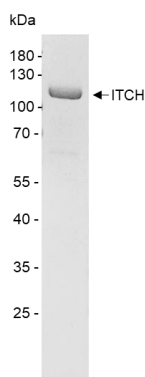
Background

ITCH is a HECT E3, which influences immunity, the skin, and plays a role in cancer by targeting various transcription factors and growth factors. ITCH is unique for its involvement in numerous cellular processes, including immune response regulation, cell cycle control, and signal transduction pathways. Dysregulation of ITCH has been linked to various diseases, making it an intriguing target for therapeutic interventions. Understanding the functions of ITCH provides valuable insights into the intricate mechanisms governing protein turnover and cellular function.

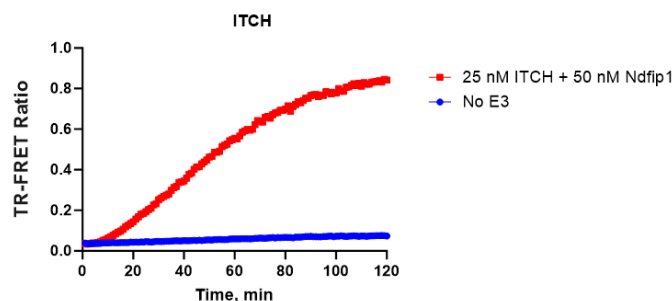
Product Information

Purity	≥ 95% by SDS-PAGE
Molecular Weight	110 kDa
Quantity	25 µg
Physical State	Liquid, 20 mM Hepes pH 7.5, 150 mM NaCl, 10% glycerol
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 ITCH. Two µg of the protein was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



Activity Assay of ITCH. 25 nM ITCH + 50 nM Ndfip1 was tested in a TR-FRET assay and shows robust signal-to-background.

References

1. Goto, J., et al., Hypertension, 2020. 76(6):1868-1878.
2. Stermer AR, et al., Experimental Biology and Medicine. 2016;241(4):367-74.

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