

Hdm2

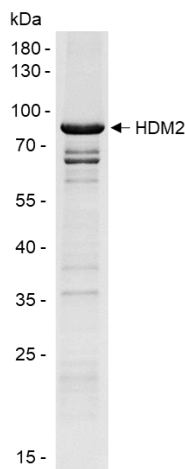
Cat. # UB313

Background

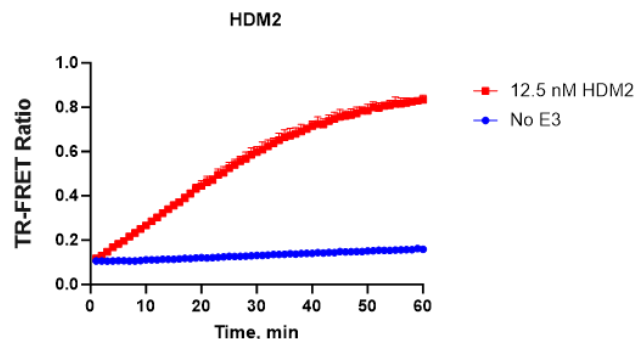
HDM2 is an E3 (ubiquitin protein ligase) that is a vital regulator of the p53 tumor suppressor protein. HDM2 can promote tumor formation by targeting tumor suppressor protein p53 for proteasomal degradation. Studies show that HDM2 has the ability to function as a ubiquitin protein ligase for both itself and p53. It has also been shown that HDM2 can mediate its E3-ligase activity in a RING finger-dependent manner. HDM2 is useful in conjunction with E1 and E2 for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Product Information

Molecular Weight	68 kDa (~80 kDa with the tag)
Quantity	20 µg
Physical State	Liquid, 1X PBS, 10% glycerol
Species	Human
Source	<i>E. coli</i>
Tag	His6-HA-Smt3
Activity	This enzyme is active with in vitro auto-ubiquitylation assay.
Storage	-80° C. Avoid repeated freeze/thaw cycles.



SDS-Page Analysis of purified HDM2. Two µg of the protein was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



Activity Assay of HDM2. 12.5 nM HDM2 was tested in a TR-FRET assay showing robust E3 ligase activity.

References

1. Koo, N., et al., Int J Mol Sci., 2022. 23(9):5005
2. Zhu, H., et al., J Hematol Oncol., 2022. 15(1):91.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2012 LifeSensors, Inc. All Rights Reserved