

Adam-10 Protease Assay Service

Adam-10 [Protease Assay](#) for compound screening and profiling via fluorescence-based quantification of enzyme activity.

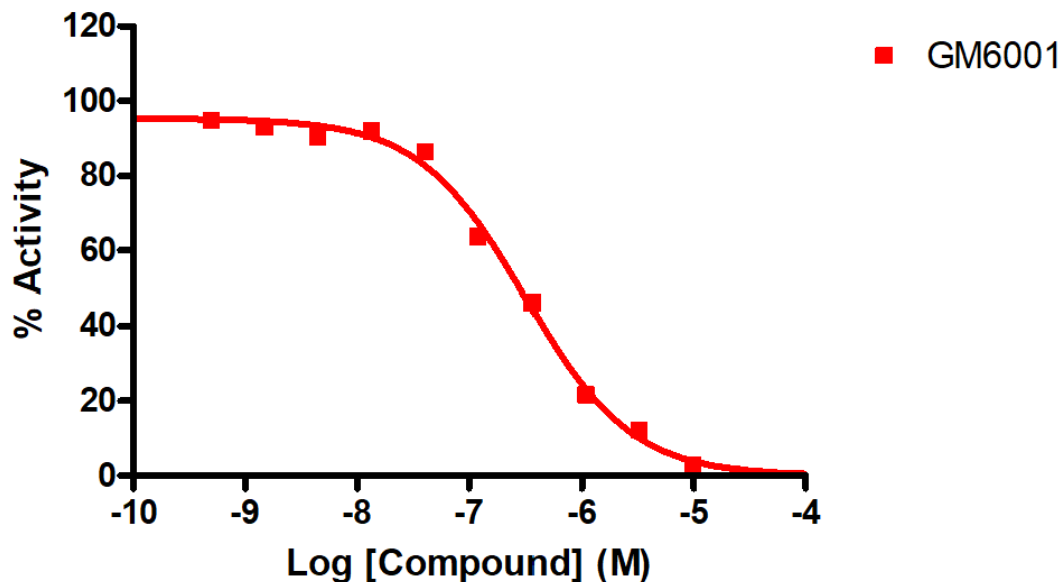
Target Overview

CAT#	Adam-10
Enzyme	Adam-10
Synonym	Disintegrin and metalloproteinase domain-containing protein 10, CDw156, Kuzbanian protein homolog, Mammalian disintegrin-metalloprotease, KUZ, MADM
Molecular Weight	52 kDa
Substrates	Fluorogenic substrate MCA-PLAQAV-Dpa-RSSSR-NH ₂ , 10 μ M

Assay Properties

Readout	Enzymatic activity
Reference Compound IC50s	Compounds IC50 (μM) GM6001 0.526
Scientific Information	O14672 (https://www.uniprot.org/uniprotkb/O14672/entry)
Screening Location	Malvern, PA, USA
Further Information	More information can be found on our website Protease Assay Services

Reference compound IC50 for Adam-10



Let's Discover Together.

© Jun 2023 **Reaction Biology**. All rights reserved.

https://www.reactionbiology.com/datasheet/adam-10_protea_malvern

