

Proteinase K Protease Assay Service

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

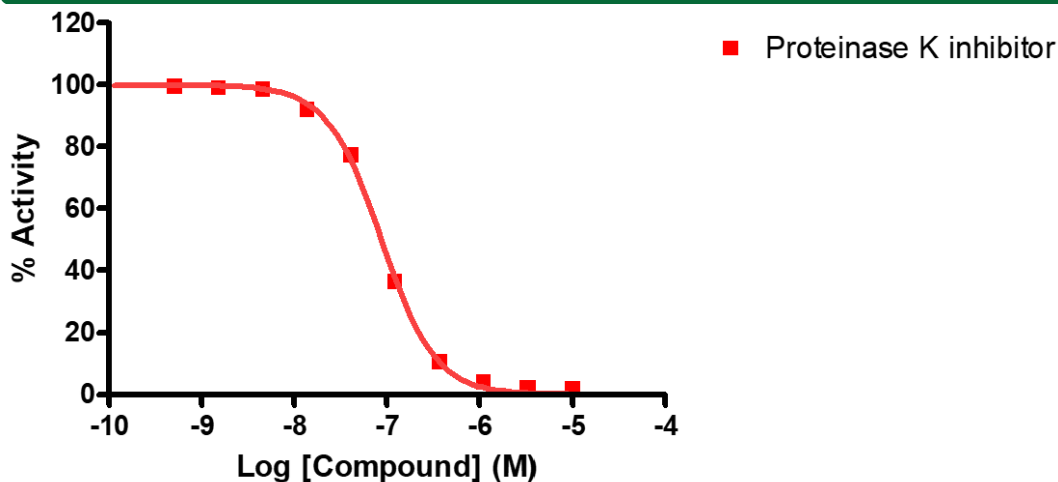
Target Overview

CAT#	Proteinase K
Enzyme	Proteinase K
Synonyms	Endopeptidase K, Tritirachium alkaline proteinase, PROK
Molecular Weight	28.9 kDa
Substrates	10 μ M fluorogenic substrate H-D-CHA-Ala-Arg-AMC.2AcOH

Assay Properties

Readout	Enzymatic activity	
Reference Compound IC50s	Compounds	IC50 (μM)
	Proteinase K Inhibitor	0.09
Scientific Information	P06873 (https://www.uniprot.org/uniprotkb/P06873/entry)	
Screening Location	Malvern, PA, USA	
Further Information	More information can be found on our website Protease Assay Services	

Reference compound IC50 for Proteinase K



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/proteinase_k_protea_malvern

