

Granzyme B Protease Assay Service

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

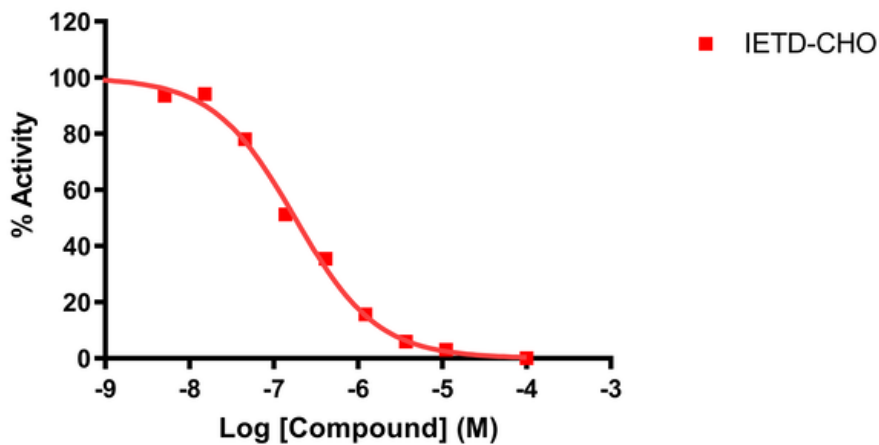
Target Overview

CAT#	Granzyme B
Enzyme	Human Granzyme B
Synonym	Granzyme-2, C11, CTLA-1, Cathepsin G-like 1, Cytotoxic T-lymphocyte proteinase 2, Fragmentin-2
Substrates	Fluorogenic substrate Z-IETD-AFC, 50 μ M

Assay Properties

Readout	Enzymatic activity
Reference Compound IC50s	Compounds IC50 (μ M) IETD-CHO 0.18
Scientific Information	P10144 (http://www.uniprot.org/uniprot/P10144)
Screening Location	Malvern, USA

Reference Compound IC50 for Granzyme B



Let's Discover Together.

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

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

