

HIV-1 Protease Assay Service

HIV-1 [Protease Assay](#) for compound screening and profiling via fluorescence-based quantification of enzyme activity.

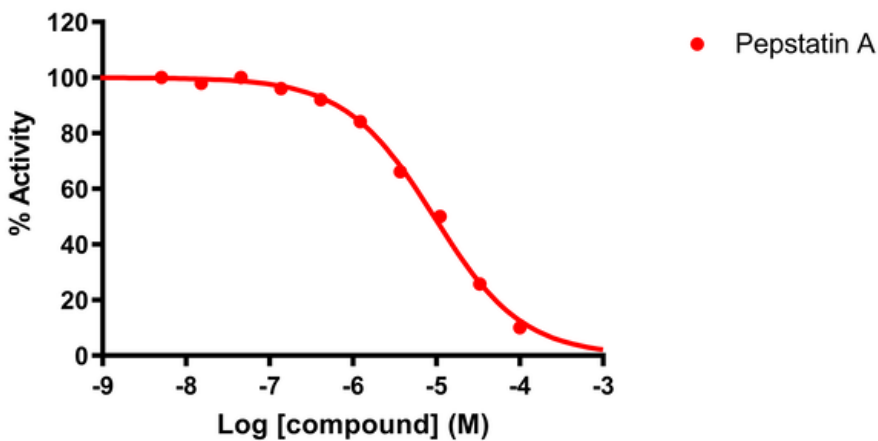
Target Overview

| | |
|------------|---|
| CAT# | HIV-1 |
| Enzyme | HIV-1 |
| Synonym | Human immunodeficiency virus-1 |
| Substrates | Fluorogenic peptide [Anaspec SensoLyte], 10 μ M |

Assay Properties

| | |
|--------------------------|---|
| Readout | Enzymatic activity |
| Reference Compound IC50s | Compounds IC50 (nM) Pepstatin A 9,100 |
| Scientific Information | NC_001802 |
| Screening Location | Malvern, USA |

Reference Compound IC50 for HIV-1



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/hiv-1_protea_malvern

