

## **Product Data Sheet**

Product Name:	HIV Protease FRET Substrate	l
Catalog Number:	AS-22992 (1 mg)	Lot Number: See label on vial
Molecular Weight: % Peak Area by HPLC:	1532.8 > 95	
Appearance:	Lyophilized red powder	
Peptide Reconstitution:	Use fresh concentrated Anhydrous DMSO as the solvent. Do not use DMSO diluted with aqueous solvents, as this is not an effective solvent. Do not add aqueous solvent to the peptide prior to DMSO addition as this may prevent proper solubilization. Add concentrated DMSO directly to the lyophilized peptide powder to obtain a final concentration of approximately 0.5mg/mL to 1mg/mL. Gently vortex to mix	
Storage:	Peptide is shipped at ambient temperature. Upon receipt, store lyophilized powder at -20°C or lower. Reconstituted peptide should be aliquoted into several freezer vials and stored at -20°C or lower. Do not freeze thaw.	
Description:	DABCYL-GABA-Ser-Gln-Asr HIV protease substrate I in s continuous assay for HIV pro- encoded by the human immu- the correct processing of vira infectious virus, and is theref acquired immunodeficiency FRET-based fluorogenic sub- site for HIV-1 PR. Incubation fluorogenic substrate resulte and a time-dependent increa related to the extent of subs- yields of the HIV- 1 PR subs- and 34.4-fold, respectively, p its simplicity and precision in for kinetic analysis, this subs- commonly used HPLC or ele substrate hydrolysis by retro Anjuere, F. et al. <i>Biochem J Lett</i> <b>262</b> , 119 (1990).	h-Tyr-Pro-IIe-Val-GIn-EDANS is also called some literature. It is widely used for the becase activity. The 11-kD protease (PR) unodeficiency virus 1 (HIV-1) is essential for al polyproteins and the maturation of fore a target for the design of selective syndrome (AIDS) therapeutics. The bestrate is derived from a natural processing n of recombinant HIV-1 PR with the ed in specific cleavage at the Tyr-Pro bond ase in fluorescence intensity that is linearly trate hydrolysis. The fluorescence quantum strate in the FRET assay increased by 40.0- per mole of substrate cleaved. Because of n the determination of reaction rates required strate offers many advantages over the ectrophoresis-based assays for peptide wiral PRs. Abs/Em = 340nm/490nm. Ref: 291, 869 (1993), Geohegan KF, et al. <i>FEBS</i>

## **Related Products:**

Name	Cat #	Size
SensoLyte® 490 HIV Protease Assay Kit *Fluorimetric*	AS-71127	1 kit
SensoLyte® 520 HIV Protease Assay Kit *Fluorimetric*	AS-71147	1 kit

## For Research Use Only