

## HIV-1 Integrase

### Human Immunodeficiency Virus 1 Integrase recombinant, *E. coli*

Cat. No.	Amount
PR-1208-1	1 mg

For *in vitro* use only  
Quality guaranteed for 12 months  
Store at -20°C

#### Avoid freeze / thaw cycles

#### Form

Liquid. Supplied in 50 mM Tris-HCl pH 8.0, 60 mM NaCl, 10 mM Glutathione, 0.25% Sarkosyl and 50% glycerol.

#### Application

May be used in ELISA and Western blots, excellent antigen for early detection of HIV seroconvertors with minimal specificity problems.

#### Specificity

Immunoreactive with all sera of HIV-1 infected individuals.

#### Molecular Weight

35 kDa including GST-Tag

#### Purity

major band and several minor smaller size bands on 12% PAGE electrophoresis followed by Coomassie staining

#### Description

The protein is a non-glycosylated polypeptide chain containing the HIV-1 immunodominant regions from the pol protein (integrase).

HIV belongs to the retrovirus family, distinguished by possession of a viral reverse transcriptase that transcribes viral RNA into DNA which is integrated into the host-cell genome.

The human immunodeficiency virus (HIV) integrase (IN), one of the three virally encoded enzymes required for HIV-1 replication, must covalently join the viral cDNA into a host chromosome for productive HIV infection.

The protein contains a GST-Tag.

#### Selected References:

Payan *et al.* (2003) [Measuring the HIV viral load with LCx (Abbott): importance for the therapeutic follow-up of 3 patients infected by type O HIV]. *Transfus. Clin. Biol.* **10**:72.

Andreola (2004) Closely related antiretroviral agents as inhibitors of two HIV-1 enzymes, ribonuclease H and integrase: "killing two birds with one stone". *Curr. Pharm. Des.* **10**:3713.

Sechi *et al.* (2004) Design and synthesis of novel indole beta-diketo acid derivatives as HIV-1 integrase inhibitors. *J. Med. Chem.* **47**:5298.

Reinke *et al.* (2004) L-chicoric acid inhibits human immunodeficiency virus type 1 integration *in vivo* and is a noncompetitive but reversible inhibitor of HIV-1 integrase *in vitro*. *Virology* **326**:203.

Hazuda *et al.* (2004) A naphthyridine carboxamide provides evidence for discordant resistance between mechanistically identical inhibitors of HIV-1 integrase. *Proc. Natl. Acad. Sci. USA* **101**:11233.

Aiello *et al.* (2004) Synthesis of novel thiazolothiazepine based HIV-1 integrase inhibitors. *Bioorg. Med. Chem.* **12**:4459.

Mousnier *et al.* (2004) Nuclear import of HIV-1 integrase is inhibited *in vitro* by styrylquinoline derivatives. *Mol. Pharmacol.* **66**:783.