





## 📕 🔳 HIV-1 p55 gag

Human Immunodeficiency Virus 1 Antigen recombinant, Baculovirus

Cat. No.	Amount
PR-1217	10 µg

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Molecular Weight: 55 kDa

Purity: > 90 % (SDS-PAGE, HPLC)

**Form:** liquid (Supplied in 10 mM Tris-HCl pH 8.0, 140 mM NaCl and 400 mM L-Arginine. For long time storage add a carrier protein like 0.1% BSA)

## **Applications:**

Antigen in ELISA and Western blots, excellent antigen for detection of HIV seroconvertors with minimal specificity problems.

## **Description:**

Recombinant HIV-1 p55 gag is glycosylated with N-linked sugars and produced in insect cells. Recombinant HIV-1 p55 is purified by proprietary chromatographic technique.

**Background:** 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. p55 is a precursor protein, which is processed to p17 (matrix), p24 (capsid), p7 (nucleocapsid), and p6 proteins by the viral protease. It is called 'assemblin' to indicate its role in viral assembly.

**Specificity:** Immunoreactive with sera of HIV-infected individuals.

## Selected References:

Ivanov *et al.* (2001) Sup35p yeast prion-like protein as an adapter for production of the Gag-p55 antigen of HIV-1 and the L-chain of botulinum neurotoxin in Saccharomyces cerevisiae. *Res. Microbiol.* **152**:27.

Neidleman *et al.* (2000) Genetically detoxified mutants of heat-labile enterotoxin from Escherichia coli are effective adjuvants for induction of cytotoxic T-cell responses against HIV-1 gag-p55. *Immunology* **101**:154.

Kazzaz *et al.* (2000) Novel anionic microparticles are a potent adjuvant for the induction of cytotoxic T lymphocytes against recombinant p55 gag from HIV-1. *J. Control. Release* **67**:347.

Bristow *et al.* (1999) Human cyclophilin has a significantly higher affinity for HIV-1 recombinant p55 than p24. *J. Acquir. Immune Defic. Syndr. Hum. Retrovirol.* **20**:334.

Jakobsen *et al.* (1998) Increased levels of soluble tumour necrosis factor receptor-I (P55) and decreased IgG1 reactivities in HIV-1 patients with cytomegalovirus disease. *Scand. J. Immunol.* **47**:591.

McAdam *et al.* (1998) Cross-clade recognition of p55 by cytotoxic T lymphocytes in HIV-1 infection. *AIDS* **12**:571.

