# **DATA SHEET**





## ■ HIV-type O Envelope

Human Immunodeficiency Virus Type O Antigen recombinant, *E. coli* 

Cat. No.	Amount
PR-1209	100 μg

For general laboratory use.

Shipping: shipped at 4 °C

Storage Conditions: store at 4 °C

Additional Storage Conditions: Please do NOT freeze

Shelf Life: 12 months

**Purity:** > 95 % (SDS-PAGE, RP-HPLC)

Form: liquid in Water

#### **Applications:**

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

#### **Description:**

Recombinant HIV type-O peptide, containing the HIV type-O transmembrane envelopederived specific sequence. Detects all clades of HIV type-O infected individuals responding to HIV type-O envelope proteins. The protein was 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. Based on genetic variability in the envelope (*env*) gene, HIV-1 can be subdivided into at least 10 distinct subtypes (designated A to J) responsible for separate geographic pandemics. Phylogenetic analyses have shown that each subtype in this major group (group M) is approximately equidistant from the others, as if arising from a common ancestor. In contrast, a few divergent HIV-1 strains form a cluster distinct from group M and have been categorized as members of the outlier group (group O).

**Specificity:** Immuno reactive with all sera of HIV type-O infected individuals.

### 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.

Miguel *et al.* (1998) Analysis of *pol* Gene Heterogeneity, Viral Quasispecies, and Drug Resistance in Individuals Infected with Group O Strains of Human Immunodeficiency Virus Type 1. *J. Virol.* **72**:9002.