

RANTES (CCL5)

Recombinant human protein expressed in E. coli cells

Catalog # R801-40N

Lot # S358-4

Product Description

Recombinant human RANTES was expressed in E. coli cells and purified by proprietary chromatographic techniques. The protein accession number is [AAH08600](#).

Gene Aliases

CCL5; SISd; SCYA5; TCP228; D17S136E; MGC17164

Endotoxin Level

Less than 0.1 ng/µg (IEU/µg) of RANTES.

Formulation

Recombinant protein stored in 1 X PBS, 10% glycerol.

Storage and Stability

Stored at desiccated below -20°C. It is stable at 4°C between 2-7 days. For long term storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

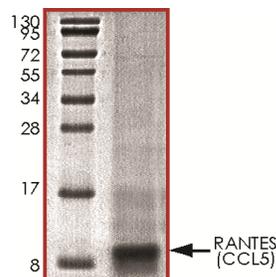
Scientific Background

RANTES is one of several CC cytokine genes clustered on the q-arm of chromosome 17 which are a family of secreted proteins involved in immunoregulatory and inflammatory processes. It causes the release of histamine from basophils and activates eosinophils. It also involved in immunomodulatory functions together with CCL5 (1). RANTES is one of the major HIV-suppressive factors produced by CD8+ cells which functions as one of the natural ligands for the chemokine receptor CCR5 and it suppresses in vitro replication of the R5 strains of HIV-1, which use CCR5 as a coreceptor (2).

References

- Hartmann, T. N. et.al: Human B cells express the orphan chemokine receptor CCR5 in a maturation-stage-dependent and CCL5-modulated manner. *Immunology* 125: 252-262, 2008.
- Arenzana-Seisdedos, F. et.al: HIV blocked by chemokine antagonist. (Letter) *Nature* 383: 400 only, 1996.

Purity



The purity was determined to be **>95%** by densitometry analysis in reducing SDS-PAGE Coomassie blue gel. Approx. MW of the major band was **8.4 kDa (7,982 Daltons)** by MW calculation).

RANTES (CCL5)

Recombinant human protein expressed in E. coli cells

Catalog Number	R801-40N
Specific Lot Number	S358-4
Purity	>95%
Concentration	0.2 µg/µl
Stability	1 yr At -20°C from date of shipment
Storage & Shipping	Stored at desiccated below -20°C. It is stable at 4°C between 2-7 days. For long term storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.