IL-8 (human),(recombinant)

Interleukin-8 (IL-8) is a chemokine, a member of the cytokine family that displays chemotactic activity for specific types of leukocytes. IL-8 regulates leukocyte recruitment and trafficking and also has mitogenic and angiogenic properties. IL-8 expression has been demonstrated in neuroendocrine tumor cells in human cancer tissue.

Ordering Information

Order Online »

ENZ-PRT185-0025

25µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Short Term Storage -20°C

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Interleukin-8, CXCL8, Monocyte-derived neutrophil

chemotactic factor, MDNCF, Neutrophil activating factor,

NAF, NAP-1

Formulation Lyophilized with no additives.

MW ~8.4 kDa

Purity ≥98% (SDS & HPLC)

Purity Detail Sterile filtered through a 0.2 micron filter.

Source Produced in *E. coli*.

UniProt ID P10145

Last modified: May 29, 2024

