IL-20 (human), (recombinant)

Interleukin 20 (IL-20) is structurally related to interleukin 10 (IL-10) and is produced by keratinocytes and monocytes. IL-20 acts through the STAT3 signaling pathway to regulate the proliferation of keratinocytes during epidermal inflammation.

Ordering Information

Order Online »

ENZ-PRT247-0010	10µg
ENZ-PRT247-0100	100μg

Manuals, SDS & CofA

View Online »

Handling & Storage

Handling Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C. For

long term storage, it is recommended to dilute to working aliquots in a 0.1% BSA

solution.

Long Term Storage -80°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Interleukin-20, Cytokine Zcyto10

Endotoxin Content ≤1 EU/µg protein (LAL test)

Formulation Lyophilized from a sterile (0.2 micron) filtered solution

containing 10 mM sodium phosphate, pH 7.5.

MW Dimer, ~17.7/35.3 kDa

Purity ≥95% (SDS-PAGE)

Reconstitution Reconstitute in sterile water at 0.1 mg/mL.

Produced in E. coli. Contains 153/306 amino acids. Source

UniProt ID Q9NYY1

Last modified: May 29, 2024

info-



eu@enzolifesciences.com