FGF-10 (human), (recombinant)

Fibroblast growth factor 10 (FGF-10) is important during embryonic development, especially during kidney, lung, limb, brain,and heart morphogenesis. FGF-10 is expressed in mesenchymal cells and facilitates epithelial-mesenchymal signaling through binding to the FGF receptor 2b (FGFR2b). FGF-10 induces the migration and invasion of cancer cells.

Ordering Information

Order Online »

ENZ-PRT234-0025	25µg
ENZ-PRT234-0100	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 Fibroblast growth factor 10, Keratinocyte growth factor 2

Biological Activity 4MBr-5 cell proliferation; ED50 \leq 200 ng/mL (\geq 5.0 x 10³

units/mg).

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 ~19.3 kDa

Purity ≥95% (SDS-PAGE)

Reconstitution Reconstitute in sterile water at 0.1 mg/mL.

Source Produced in *E. coli*. Contains 170 amino acids.

UniProt ID O15520

Last modified: May 29, 2024

