

DESCRIPTION

Source *E. coli*-derived FGF-10 protein
Accession # O15520.1

Predicted Molecular Mass 17 kDa

SPECIFICATIONS

SDS-PAGE Monomeric FGF-10 protein only

Activity No significant difference between EC₅₀ of reference and test lots

Endotoxin Level <0.10 EU per 1 µg of the protein by the LAL method.

Mass Spectrometry Single species with expected mass

Mycoplasma Negative when tested in both ribosomal RNA hybridization and luminescence assays

Formulation Lyophilized from HEPES/NaCl/mannitol See Certificate of Analysis for details.

PREPARATION AND STORAGE

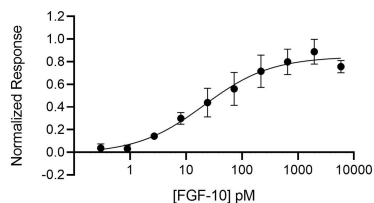
Reconstitution Resuspend in water at >100 µg/ml, prepare single use aliquots, add carrier protein if desired.

Shipping The product is shipped lyophilized at ambient temperature, on ice blocks or on dryice. Shipping at ambient temperature does not affect the bioactivity or stability of the protein. Upon receipt, store immediately at the conditions stated below.

Stability & Storage BulkLotPrefix assignment required for Storage Info

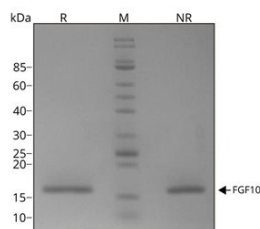
DATA

Bioactivity



Recombinant Human/Rat/Bovine FGF-10, Animal-Free Protein Bioactivity
FGF-10 activity is determined using the firefly luciferase reporter assay (*) in stably transfected HEK293T cells. Cells are treated in triplicate with a serial dilution of FGF-10. Firefly luciferase activity is measured and normalized. EC₅₀ = 0.36 ng/ml (21.1 pM).
*Promega pGL4.33[luc2P/SRE/Hygro] #E1340

N/A



Recombinant Human/Rat/Bovine FGF-10, Animal-Free Protein SDS-PAGE
FGF-10 migrates as a single band at 17 kDa in non-reducing (NR) conditions and upon reduction (R). No contaminating protein bands are visible. Purified recombinant human FGF-10 protein (7 µg) was resolved using 15% w/v SDS-PAGE in reduced (+β-mercaptoethanol, R) and non-reduced conditions (NR) and stained with Coomassie Brilliant Blue R250.

BACKGROUND

The Fibroblast Growth Factors (FGFs) are heparin binding glycoproteins that exert a variety of biological activities toward cells of mesenchymal, neuronal, and epithelial origin. FGF-10 belongs to the subgroup of FGFs that also includes FGF-3, -7, and -22 (1). Mature human FGF-10 is an approximately 20 kDa protein that contains a serine-rich region near its N-terminus (2, 3). It shares 93% and 96% amino acid sequence identity with mouse and rat FGF-10, respectively. FGF-10 is secreted by mesenchymal cells and associates with extracellular FGF-BP (1, 4). It preferentially binds and activates epithelial cell FGF R2 (IIIb) and interacts more weakly with FGF R1 (IIIb) (5). The mitogenic and chemotactic properties of FGF-10 are critical in many tissues during embryogenesis. This includes limb bud initiation (6), palate development (7), branching morphogenesis and directional outgrowth of lung buds (8, 9), formation of the otic vesicle and chochlea (10), adipogenesis (11), and the development of prostate, mammary, lacrimal, and submandibular salivary glands (12 - 15). FGF R2 (IIIb) signaling in these responsive tissues is similarly important during embryogenesis (7, 10, 13 - 15). The expression and function of FGF-10 are negatively regulated by Shh and BMP-4 in the developing lung (8, 9). Overlapping expression patterns and activities with FGF-3, -7, and -8 suggest at least a partial redundancy in FGF-10 biology (7, 10, 14, 15). FGF-10 induced signaling through FGF R2 (IIIb) also contributes to the progression of pancreatic cancer (16).

References:

1. Beenken, A. and M. Mohammadi (2009) *Nat. Rev. Drug Discov.* **8**:235.
2. Igarashi, M. *et al.* (1998) *J. Biol. Chem.* **273**:13230.
3. Emoto, H. *et al.* (1997) *J. Biol. Chem.* **272**:23191.
4. Beer, H.-D. *et al.* (2005) *Oncogene* **24**:5269.
5. Zhang, X. *et al.* (2006) *J. Biol. Chem.* **281**:15694.
6. Min, H. *et al.* (1998) *Genes Dev.* **12**:3156.
7. Rice, R. *et al.* (2004) *J. Clin. Invest.* **113**:1692.
8. Bellusci, S. *et al.* (1997) *Development* **124**:4867.
9. Weaver, M. *et al.* (2000) *Development* **127**:2695.
10. Pirvola, U. *et al.* (2000) *J. Neurosci.* **20**:6125.
11. Sakaue, H. *et al.* (2002) *Genes Dev.* **16**:908.
12. Donjacour, A.A. *et al.* (2003) *Dev. Biol.* **261**:39.
13. Mailloux, A.A. *et al.* (2002) *Development* **129**:53.
14. Makarenkova, H.P. *et al.* (2000) *Development* **127**:2563.
15. Jaskoll, T. *et al.* (2005) *BMC Dev. Biol.* **5**:11.
16. Nomura, S. *et al.* (2008) *Br. J. Cancer* **99**:305.

PRODUCT SPECIFIC NOTICES

The above product was manufactured, tested and released by R&D System's contract manufacturer, Qkine Ltd, at 1 Murdoch House, Cambridge, UK, CB5 8HW. The product is for research use only and not for the diagnostic or therapeutic use.