

PRODUCT INFORMATION



FGF7 (human, recombinant)

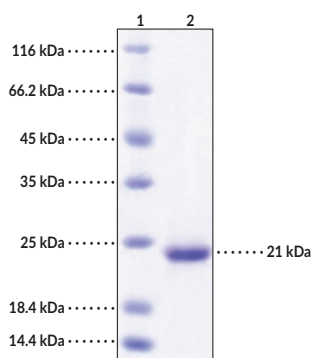
Item No. 42756

Overview and Properties

Synonyms:	Fibroblast Growth Factor 7, HBGF-7, Heparin-binding Growth Factor 7, Keratinocyte Growth Factor
Source:	Active recombinant human N-terminal His-tagged FGF7 expressed in <i>E. coli</i>
Amino Acids:	32-194
Uniprot No.:	P21781
Molecular Weight:	21 kDa
Storage:	-80°C (as supplied)
Stability:	≥1 year
Purity:	≥95% estimated by SDS-PAGE
Supplied in:	Lyophilized from sterile 50 mM Tris, pH 7.5, with 250 mM sodium chloride, 5% trehalose, 5% mannitol, and 0.01% polysorbate 80
Endotoxin Testing:	< 1.0 EU/μg, determined by the LAL endotoxin assay

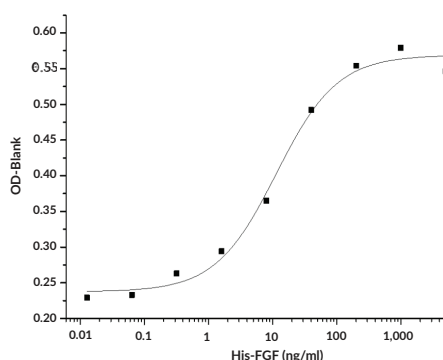
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



Lane 1: MW Markers
Lane 2: FGF7

SDS-PAGE Analysis of FGF7. This protein has a calculated molecular weight of 21 kDa.



Cell proliferation assay using using BaF3 mouse pro-B cells transfected with human FGFR2b. The EC₅₀ value for this effect is typically 3-20 ng/ml.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/01/2026

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Fibroblast growth factor 7 (FGF7), also known as keratinocyte growth factor (KGF), is a member of the FGF7 subfamily.^{1,2} It is composed of an N-terminal signaling region, which is cleaved prior to secretion, a Ca backbone, a heparin-binding domain, and a receptor-binding domain.^{2,3} FGF7 is secreted by stromal fibroblasts and forms a quaternary complex with FGFR2b dimers, and to a lesser extent FGFR1b dimers, with heparan sulfate proteoglycans to induce ERK- or MAPK-induced proliferation in epithelial cells.^{1,3,4} It is also necessary for proper nephron and hippocampal development and post-radiation thymic regeneration in mice.⁴ Knockout of *Fgf7* decreases cartilage erosion, osteophyte formation, bone density, and subchondral bone plate thickness in a mouse model of osteoarthritis induced by destabilization of the medial meniscus (DMM).⁵ Increased levels of FGF7 in cancer-associated fibroblasts are correlated with increased arterial and lymphatic metastasis, increased amount of residual tumor following surgery, and advanced cancer staging in patients with ovarian cancer.⁶ Cayman's FGF7 (human, recombinant) protein can be used for cell-based assays. This protein consists of 162 amino acids and has a calculated molecular weight of 21 kDa.

References

1. Mason, I. Initiation to end point: The multiple roles of fibroblast growth factors in neural development. *Nat. Rev. Neurosci.* **8**(8), 583-596 (2007).
2. Ye, S., Luo, Y., Lu, W., *et al.* Structural basis for interaction of FGF-1, FGF-2, and FGF-7 with different heparan sulfate motifs. *Biochemistry* **40**(48), 14429-14439 (2001).
3. Finch, P.W., Rubin, J.S., Miki, T., *et al.* Human KGF is FGF-related with properties of a paracrine effector of epithelial cell growth. *Science* **245**(4919), 752-755 (1989).
4. Zinkle, A. and Mohammadi, M. Structural biology of the FGF7 subfamily. *Front. Genet.* **10**, 102 (2019).
5. Shin, Y., Kwak, J.-S., Kim, S.K., *et al.* Fibroblast growth factor 7 (FGF7) causes cartilage destruction, subchondral bone remodeling, and the premature growth plate closure in mice. *Osteoarthritis Cartilage* (2024).
6. Feng, S., Ding, B., Dai, Z., *et al.* Cancer-associated fibroblast-secreted FGF7 as an ovarian cancer progression promoter. *J. Transl. Med.* **22**(1), 280 (2024).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM