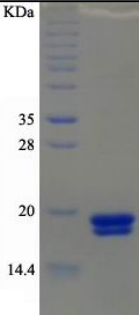


Recombinant Human FGF2 Protein Data Sheet

Catalog #	hRP-L0207-EF011
Size	25 µg
Protein Name	Human fibroblast growth factor 2
Protein Symbol	FGF2
Original Source	<i>Homo sapiens</i>
Expression System	<i>E.coli</i>
GenBank Accession #	M27968.1
Uniprot Accession #	P09038-2
Description	FGF2 which is a member of the fibroblast growth factor (FGF) family. FGF family members bind heparin and possess broad mitogenic and angiogenic activities. This protein has been implicated in diverse biological processes, such as limb and nervous system development, wound healing, and tumor growth.
Application	WB, ELISA, IP, antibody production, protein array
Fusion tag	N-His
Peptide Length	170aa(including fusion tag)
Molecular Weight	19.2kDa(including fusion tag)
pI	9.9
Activity	NA
Storage	Storage buffer: 50mM Tris.Cl, 50mM NaCl, 2.5% Glycerol, pH8.0. Store at -80°C and avoid repeated freeze-thaw cycles.
	
Reference:	• Memory of fearful events: the role of fibroblast growth factor-2 in fear acquisition and extinction • The discovery of basic fibroblast growth factor/fibroblast growth factor-2 and its role in haematological malignancies • Fibroblast growth factor-2 and cardioprotection • A local application of recombinant human fibroblast growth factor 2 for tibial shaft fractures: A randomized, placebo-controlled trial • Fibroblast growth factor 2: from laboratory evidence to clinical application

GeneCopoeia™
Expressway to Discovery

GeneCopoeia Inc.

9620 Medical Center Drive, Suite 101

Rockville, MD 20850

USA

Phone: 301-762-0888

Toll free: 1-301-762-0888

Fax: 301-762-3888

Web: www.genecopoeia.com

Inquiry: inquiry@genecopoeia.com

Technical Support: support@genecopoeia.com