

Basic Fibroblast Growth Factor (human) (recombinant)

Ordering Information

[Order Online »](#)

RLT-300-003	50µg
-------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	After reconstitution, prepare aliquots and store at -20°C. Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	bFGF (human) (recombinant), FGF-2 (human) (recombinant)
Biological Activity	Stimulates ³ H-thymidine incorporation in primary human macro- or microvascular endothelial cells (ED ₅₀ =0.4-1.0ng/ml). Maximal effect for DNA synthesis was observed at 1.0-2.0ng/ml.
Endotoxin Content	≤1.2EU/μg protein.
Formulation	Lyophilized.
MW	~16kDa.
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute with 50mM acetic acid, containing at least 0.1% HSA or BSA (>10μg/ml).
Source	Produced in <i>E. coli</i> .
Specific Activity	~2.5×10 ⁶ U/mg.
Technical Info / Product Notes	For the Original Manufacturer's data sheet please click here .



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)