

Vascular Endothelial Growth Factor C (C152S mutant) (rat) (recombinant) (His)

This mutant is lacking binding to VEGFR-2 (KDR) and can serve as a unique tool to investigate the signal transduction of VEGF-C via VEGFR-3 (Flt-4). According to the high degree of homology between growth factors of mammalian species, the protein can also be used for experiments with human cells and for further *in vitro* and *in vivo* experiments with other cells and species.

Ordering Information

[Order Online »](#)

RLT-R20-017	20µg
-------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	After reconstitution, prepare aliquots and store at -20°C. Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	VEGF-C (C152S mutant) (rat) (recombinant) (His)
Biological Activity	Stimulates phosphorylation of VEGFR-3 in pig aortic endothelial cells (ED ₅₀ =150-300ng/ml).
Endotoxin Content	≤1.2EU/μg protein.
Formulation	Lyophilized.
MW	~18-24kDa.
Purity	≥90% (SDS-PAGE)
Reconstitution	Reconstitute with PBS or medium (>50μg/ml).
Solubility	Soluble in water and most aqueous buffers.
Source	Produced in insect cells.
Specific Activity	~5×10 ³ 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

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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