

Growth Hormone, placental 22K (human), (recombinant)

Ordering Information

[Order Online »](#)

RLT-500-007	50µg
-------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Growth hormone variant, GH-V, Growth hormone 2, Placenta-specific growth hormone
Biological Activity	rHuGH-22K-pl is fully biologically active when compared to WHO reference standard using in vitro bioassay. It is also capable of forming a 1:2 complex with the recombinant human growth hormone receptor extracellular domain (ECD).
Endotoxin Content	< 0.1 ng/μg of protein (<1EU/μg)
Formulation	Lyophilized.
MW	22.3 kDa
Purity	> 98.0% as determined by RP-HPLC, Gel filtration and SDS-PAGE Silver Stained gel.
Source	Produced in <i>E. coli</i> .
UniProt ID	P01242-2



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