

Connective tissue growth factor (human), (recombinant)

Connective Tissue Growth Factor (CTGF) is a member of cysteine rich regulatory proteins that are both mitogenic and chemotactive. CTGF has numerous effects on development and differentiation.

Ordering Information

[Order Online »](#)

ENZ-PRT212-0020	20µg
ENZ-PRT212-0100	100µg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C. For long term storage, it is recommended to dilute to working aliquots in a 0.1% BSA solution.
Short Term Storage	+4°C
Long Term Storage	-80°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	CCN2, HCS24, CTGF
Endotoxin Content	≤0.1 EU/μg using USP 85 / EP 2.6.14
Formulation	Lyophilized from a 0.2 μm filtered solution containing 0.1% Trifluoroacetic Acid (TFA).
MW	~11.2 kDa
Purity	≥90% using Coomassie Staining
Reconstitution	Reconstitute in sterile 5 mM sodium acetate, pH 6.0, at 0.1 mg/mL.
Source	Produced in <i>E. coli</i> . A non-glycosylated protein containing 98 amino acids.
UniProt ID	P29279

Last modified: June 17, 2025



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-
fr@enzolifesciences.com](mailto:info-fr@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)