

SARTORIUS

Stock Up and Save on Research Grade Growth Factors & Cytokines

Save 30%!



Sartorius' range of high-quality research use only (RUO) growth factors and cytokines are manufactured to the highest quality standards using recombinant DNA technology which do not contain any animal derived components or contaminants.

Stock up and save! Get a 30% discount on RUO cytokines and growth factors now through May 31, 2026. Applicable in eligible countries.

Just enter the code **CYTOK2026** on the eShop to take advantage of this offer.

- Promotion valid until May 31, 2026.
- Applicable in EMEA and NA
- Cannot be combined with any other offers or discounts



Research Grade Growth Factors & Cytokines Features

SKU Number—Material Description

CYK-0010-0003 - Recomb human KGF (FGF-7) protein 0.01mg
CYK-0010-0005 - Recomb human VEGF 165 protein 0.01mg
CYK-0010-0008 - Recombinant human BDNF protein 0.01mg
CYK-0025-0003 - Recomb human KGF (FGF-7) protein 0.025mg
CYK-0025-0005 - Recomb human VEGF 165 protein 0.025mg
CYK-0025-0006 - Recombinant human BMP-4 protein 0.025mg
CYK-0025-0007 - Recomb human Activin A protein 0.025mg
CYK-0025-0008 - Recombinant human BDNF protein 0.025mg
CYK-0025-0009 - Recombinant human GDNF protein 0.025mg
CYK-0025-0010 - Recombinant human TGF-β1 PLUS protein
CYK-0025-0011 - Recomb human PDGF-AA protein 0.025mg
CYK-0025-0012 - Recombinant human LIF protein 0.025mg
CYK-0025-0013 - Recombinant human BMP-2 protein 0.025mg
CYK-0025-0014 - Recombinant human NRG-1 protein 0.025mg
CYK-0025-0015 - Recombinant human TGF-β3 PLUS protein
CYK-0025-0016 - Recombinant human Noggin protein 0.025mg
CYK-0025-0017 - Recombinant human FGF-10 protein 0.025mg
CYK-0025-0018 - Recombinant human FGF-9 protein 0.025mg
CYK-0025-0019 - Recomb human R-spondin 1 protein 0.025mg
CYK-0025-0020 - Recombinant human NT-3 protein 0.025mg
CYK-0100-0021 - Recomb human HGF (NK1) protein 0.1mg
CYK-0025-0022 - Recombinant human OSM protein 0.025mg
CYK-0025-0024 - Recombinant human SHH protein 0.025mg
CYK-0025-0025 - Recombinant human FGF-4 protein 0.025mg
CYK-0025-0026 - Recomb human R-spondin 3 protein 0.025mg
CYK-0025-0027 - Recombinant human FGF-8a protein 0.025mg
CYK-0025-0028 - Recombinant human FGF-8b protein 0.025mg
CYK-0025-0030 - Recombinant human G-CSF protein 0.025mg
CYK-0025-0031 - Recombinant human M-CSF protein 0.025mg
CYK-0025-0032 - Recomb human PDGF-BB protein 0.025mg
CYK-0025-1001 - Recombinant Human Interleukin-2
CYK-0025-1002 - Recombinant Human Stem Cell Factor
CYK-0025-1003 - Recombinant Human Interleukin-15
CYK-0025-1004 - Recombinant human Tumour Necrosis Factor
CYK-0025-1005 - Recombinant Human Fms-like tyrosine kina
CYK-0025-1006 - Recombinant Human Interleukin-4
CYK-0025-1007 - Recombinant Human Interleukin-10
CYK-0025-1008 - Recombinant Human Interleukin-7
CYK-0025-1009 - Recombinant Human Interleukin-12
CYK-0025-1011 - Recombinant Human Interleukin-3
CYK-0025-1012 - Recombinant Human Thrombopoietin
CYK-0025-1013 - Recombinant Human Erythropoietin
CYK-0025-1014 - Recombinant Human Interleukin-6
CYK-0025-1015 - Recombinant Human Interleukin-18
CYK-0025-1016 - Recombinant Human Interleukin-1 Beta
CYK-0025-1017 - Recombinant Human Interferon gamma
CYK-0025-1018 - Recombinant Human Interleukin-21
CYK-0025-1019 - Recombinant Human Granulocyte-macrophage
CYK-0050-0002 - Recombinant FGF2-G3 protein 0.05mg
CYK-0050-0003 - Recombinant human KGF (FGF-7) protein
CYK-0050-0010 - Recombinant human TGF-β1 PLUS protein
CYK-0050-0015 - Recombinant human TGF-β3 PLUS protein
CYK-0050-0020 - Recombinant human NT-3 protein 0.05mg
CYK-0100-0001 - Recombinant human EGF protein 0.1mg
CYK-0100-0002 - Recombinant FGF2-G3 protein 0.1mg
CYK-0100-0004 - Recomb human IGF-1 LR3 protein 0.1mg
CYK-0100-0005 - Recomb human VEGF 165 protein 0.1mg

SKU Number—Material Description

CYK-0100-0006 - Recombinant human BMP-4 protein 0.1mg
CYK-0100-0007 - Recomb human Activin A protein 0.1mg
CYK-0100-0008 - Recombinant human BDNF protein 0.1mg
CYK-0100-0009 - Recombinant human GDNF protein 0.1mg
CYK-0100-0010 - Recombinant human TGF-β1 PLUS protein
CYK-0100-0011 - Recomb human PDGF-AA protein 0.1mg
CYK-0100-0012 - Recombinant human LIF protein 0.1mg
CYK-0100-0013 - Recombinant human BMP-2 protein 0.1mg
CYK-0100-0014 - Recombinant human NRG-1 protein 0.1mg
CYK-0100-0015 - Recombinant human TGF-β3 PLUS protein
CYK-0100-0016 - Recombinant human Noggin protein 0.1mg
CYK-0100-0017 - Recombinant human FGF-10 protein 0.1mg
CYK-0100-0018 - Recombinant human FGF-9 protein 0.1mg
CYK-0100-0019 - Recomb human R-spondin 1 protein 0.1mg
CYK-0100-0020 - Recombinant human NT-3 protein 0.1mg
CYK-0025-0021 - Recomb human HGF (NK1) protein 0.025mg
CYK-0100-0022 - Recombinant human OSM protein 0.1mg
CYK-0100-0023 - Recomb human FGF-2 protein 154 aa 0.1mg
CYK-0100-0024 - Recombinant human SHH protein 0.1mg
CYK-0100-0025 - Recombinant human FGF-4 protein 0.1mg
CYK-0100-0026 - Recomb human R-spondin 3 protein 0.1mg
CYK-0100-0027 - Recombinant human FGF-8a protein 0.1mg
CYK-0100-0028 - Recombinant human FGF-8b protein 0.1mg
CYK-0100-0030 - Recombinant human G-CSF protein 0.1mg
CYK-0100-0031 - Recombinant human M-CSF protein 0.1mg
CYK-0100-0032 - Recomb human PDGF-BB protein 0.1mg
CYK-0100-1001 - Recombinant Human Interleukin-2
CYK-0100-1002 - Recombinant Human Stem Cell Factor
CYK-0100-1003 - Recombinant Human Interleukin-15
CYK-0100-1004 - Recombinant human Tumour Necrosis Factor
CYK-0100-1005 - Recombinant Human Fms-like tyrosine kina
CYK-0100-1006 - Recombinant Human Interleukin-4
CYK-0100-1007 - Recombinant Human Interleukin-10
CYK-0100-1008 - Recombinant Human Interleukin-7
CYK-0100-1009 - Recombinant Human Interleukin-12
CYK-0100-1011 - Recombinant Human Interleukin-3
CYK-0100-1012 - Recombinant Human Thrombopoietin
CYK-0100-1013 - Recombinant Human Erythropoietin
CYK-0100-1014 - Recombinant Human Interleukin-6
CYK-0100-1015 - Recombinant Human Interleukin-18
CYK-0100-1016 - Recombinant Human Interleukin-1 Beta
CYK-0100-1017 - Recombinant Human Interferon gamma
CYK-0100-1018 - Recombinant Human Interleukin-21
CYK-0100-1019 - Recombinant Human Granulocyte-macrophage
CYK-0500-0001 - Recombinant human EGF protein 0.5mg
CYK-0500-0002 - Recombinant FGF2-G3 protein 0.5mg
CYK-0500-0003 - Recomb human KGF (FGF-7) protein 0.1mg
CYK-0500-0004 - Recomb human IGF-1 LR3 protein 0.5mg
CYK-0500-0005 - Recomb human VEGF 165 protein 0.5mg
CYK-0500-0007 - Recomb human Activin A protein 0.5mg
CYK-0500-0008 - Recombinant human BDNF protein 0.5mg
CYK-0500-0009 - Recombinant human GDNF protein 0.5mg
CYK-0500-0014 - Recombinant human NRG-1 protein 0.5mg
CYK-0500-0016 - Recombinant human Noggin protein 0.5mg
CYK-0500-0017 - Recombinant human FGF-10 protein 0.5mg
CYK-0500-0019 - Recomb human R-spondin 1 protein 0.5mg
CYK-0500-0023 - Recomb human FGF-2 protein 154 aa 0.5mg

Terms & Conditions:

- Discount applies to current list price specified for your country, excluding sales tax.
- Cannot be combined with other offers or discounts. Valid until May 31, 2026.

- Promotion code must be provided at time of order.
- **Eligible Countries:** Austria, Belgium, Canada, Denmark, Finland, France, Germany, Iceland, Ireland, Italy, Netherlands, Norway, North America, Portugal, Spain, Sweden, Switzerland, United Kingdom

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
3874 Research Park Drive
Ann Arbor, MI 48108
Phone +1 734 769 1600