

Exodus-2 (human), (recombinant)

Exodus-2 (CCL21 or 6CKine), is a chemokine produced in the lymph nodes and spleen. Exodus-2 signals through the chemokine receptor CCR7 to regulate thymocyte and activated T cell migration and also mediates the homing of lymphocytes to the lymphatic system. Human and mouse Exodus-2 share > 85% amino acid identity.

Ordering Information

[Order Online »](#)

ENZ-PRT198-0020	20µg
ENZ-PRT198-0100	100µg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Reconstituted solution is stable for 3 months when stored at -20°C to -80°C and for 1 month when stored at 4°C after reconstituting as directed.
Handling	Centrifuge vial before opening. Suspend the product by gently pipetting the above recommended solution and allow several minutes for complete reconstitution. Do not vortex. For prolonged storage, dilute to working aliquots in a 0.1% BSA solution, store at -80°C. Avoid freeze/thaw cycles.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	CCL21, 6Ckine
Endotoxin Content	≤1 EU/μg (LAL)
Formulation	Lyophilized from a sterile (0.2 micron) filtered aqueous solution containing 0.1% trifluoroacetic acid (TFA).
MW	Monomer, 12.3 kDa (111 amino acids)
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute with sterile water at 0.1 mg/mL.
Sequence	SDGGAQD CCLKYSQRKI PAKVVRSYRK QEPSLGCSIP AILFLPRKRS QAELCADPKE LWVQQLMQHL DKTPSPQKPA QGCRKDRGAS KTGKKKGKGSK GCKRTERSQT PKG
Source	Produced in <i>E. coli</i> .
UniProt ID	O00585



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)