# 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 KTGKKGKGSK

**GCKRTERSQT PKG** 

Source Produced in *E. coli*.



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com O00585
Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

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