

ZK-756326

CCR8 activator

Novel small molecule agonist at the CCR8 chemokine receptor (EC_{50} =254nM) with no activity at CCR1, CCR2B, CCR4, CCR5, CXCR2, CXCR3 or CXCR4. Dose dependently increases intracellular calcium and stimulates extracellular acidification in cells expressing human CCR8. Murine CCR8-expressing cells underwent chemotaxis on exposure to this agonist. Inhibits HIV fusion of cells expressing CD4 and CCR8. Small molecule CCR8 agonists represent useful tools for exploring CCR8 biology especially in experiments where the use of the natural protein agonist is impractical.

Citations: 2

[View Online »](#)

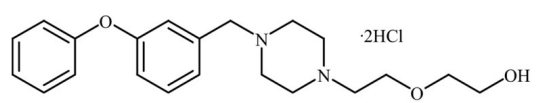
Ordering Information

[Order Online »](#)

BML-CX101-0010	10mg
BML-CX101-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at room temperature for up to one year. Store solutions at -20°C for up to 3 months.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2-[2-[4-[(3-Phenoxyphenyl)methyl]-1-piperazinyl]ethoxy]ethanol dihydrochloride
Appearance	White solid.
CAS	874911-96-3
Couple Target	Chemokine receptor
Couple Type	Activator, Ligand
Formula	$C_{21}H_{28}N_2O_3 \cdot 2HCl$
MW	356.4 . 73.0
Purity	≥98% (TLC)
Solubility	Soluble in water (5 mg/ml) or DMSO (4 mg/ml).



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-fr@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