

# 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 »](#)

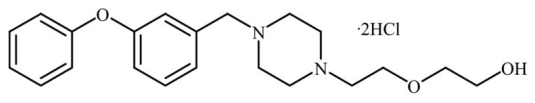
## Ordering Information

[Order Online »](#)

BML-CX101-0010	10mg
----------------	------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	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.
<b>Long Term Storage</b>	Ambient
<b>Shipping</b>	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

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

Last modified: May 29, 2024



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