

CITCO

Constitutive androstane receptor ligand

CITCO is a novel potent and selective agonist for the human orphan nuclear receptor CAR (constitutive androstane receptor). CITCO displayed an EC_{50} of 49 nM in a CAR FRET assay and showed 50-fold selectivity for CAR over PXR. CITCO displayed no activity against a panel of 15 other nuclear receptors at a concentration of 10 μ M. In hepatocytes CITCO induced CAR nuclear translocation and induced the CAR target gene CYP2B6.

Citations: 6

[View Online »](#)

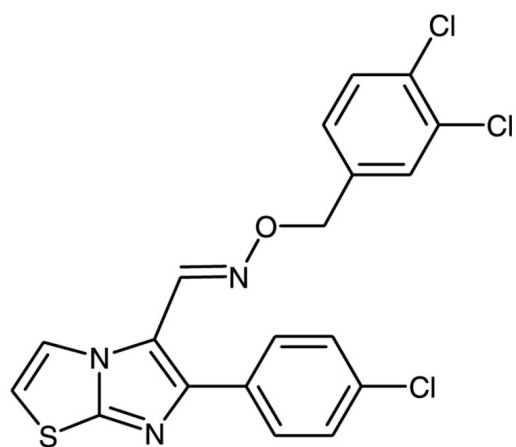
Ordering Information

[Order Online »](#)

BML-GR233-0025	25mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 6-(4-Chlorophenyl)imidazo[2,1-b] [1,3]thiazole-5-carbaldehyde O-3,4-dichlorobenzyl) oxime

Appearance Pale orange solid.

CAS 338404-52-7

Couple Target Constitutive androstane receptor

Couple Type Ligand

Formula $C_{19}H_{12}Cl_2N_3OS$

MW 436.7

Solubility Soluble in DMF (25mg/ml) or DMSO (10mg/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