

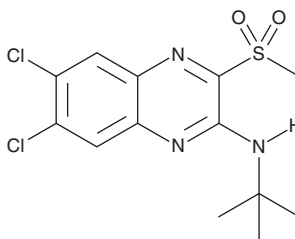
PRODUCT INFORMATION



GLP-1R Agonist DMB

Item No. 29678

CAS Registry No.: 281209-71-0
Formal Name: 6,7-dichloro-N-(1,1-dimethylethyl)-3-(methylsulfonyl)-2-quinoxalinamine
MF: C₁₃H₁₅Cl₂N₃O₂S
FW: 348.2
Purity: ≥98%
UV/Vis.: λ_{max}: 223, 264 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

GLP-1R agonist DMB is supplied as a crystalline solid. A stock solution may be made by dissolving the GLP-1R agonist DMB in the solvent of choice, which should be purged with an inert gas. GLP-1R agonist DMB is soluble in DMSO.

Description

GLP-1R agonist DMB is an agonist of glucagon-like peptide 1 receptor (GLP-1R; K_B = 26.3 nM for the recombinant human receptor).¹ It inhibits forskolin-induced cAMP accumulation in BHK cells expressing the recombinant human receptor with an EC₅₀ value of 101 nM.² GLP-1R (100 and 1,000 nM) increases insulin levels induced by D-(+)-glucose (Item No. 23733) in pancreatic islets isolated from wild-type, but not *Glp1r* knockout, mice. It reduces fasting plasma glucose levels in mice when administered at a dose of 5 μmol/kg.³

References

1. Koole, C., Wootten, D., Simms, J., *et al.* Allosteric ligands of the glucagon-like peptide 1 receptor (GLP-1R) differentially modulate endogenous and exogenous peptide responses in a pathway-selective manner: Implications for drug screening. *Mol. Pharmacol.* **78**(3), 456-465 (2010).
2. Knudsen, L.B., Kiel, D., Teng, M., *et al.* Small-molecule agonists for the glucagon-like peptide 1 receptor. *Proc. Natl. Acad. Sci. USA* **104**(3), 937-942 (2007).
3. Irwin, N., Flatt, P.R., Patterson, S., *et al.* Insulin-releasing and metabolic effects of small molecule GLP-1 receptor agonist 6,7-dichloro-2-methylsulfonyl-3-N-*tert*-butylaminoquinoxaline. *Eur. J. Pharmacol.* **628**(1-3), 268-273 (2010).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/15/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM