

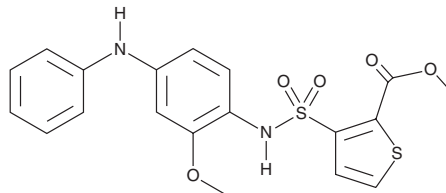
PRODUCT INFORMATION



GSK0660

Item No. 15272

CAS Registry No.: 1014691-61-2
Formal Name: 3-[[[2-methoxy-4-(phenylamino)phenyl]amino]sulfonyl]-2-thiophenecarboxylic acid, methyl ester
MF: C₁₉H₁₈N₂O₅S₂
FW: 418.5
Purity: ≥95%
UV/Vis.: λ_{max}: 286 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

GSK0660 is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK0660 in the solvent of choice. GSK0660 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of GSK0660 in these solvents is approximately 2, 12, and 20 mg/ml, respectively.

GSK0660 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GSK0660 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. GSK0660 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

GSK0660 is the first reported antagonist of PPARβ/δ (IC₅₀ = 155 nM) with no measurable affinity for PPARα or PPARγ (IC₅₀s ≥ 10 μM).¹ Because of the limited bioavailability of GSK0660, the related compound GSK3787 (Item No. 15219) was developed as a selective PPARβ/δ antagonist with more suitable pharmacokinetic properties.²

References

1. Shearer, B.G., Steger, D.J., Way, J.M., *et al.* Identification and characterization of a selective peroxisome proliferator-activated receptor β/δ (NR1C2) antagonist. *Molecular Endocrinology* **22**(2), 523-529 (2008).
2. Palkar, P.S., Borland, M.G., Naruhn, S., *et al.* Cellular and pharmacological selectivity of the peroxisome proliferator-activated receptor-β/δ antagonist GSK3787. *Mol. Pharmacol.* **78**(3), 419-430 (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, 11/16/2022

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