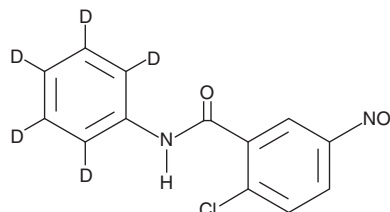


PRODUCT INFORMATION



GW 9662-d₅
Item No. 9000497

Formal Name: 2-chloro-5-nitrobenzanilide-
10,11,12,13,14-d₅
MF: C₁₃H₄ClD₅N₂O₃
FW: 281.7
Chemical Purity: ≥98% (GW 9662)
Deuterium
Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
UV/Vis.: λ_{max}: 260 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

GW 9662-d₅ is intended for use as an internal standard for the quantification of GW 9662 by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

GW 9662-d₅ is supplied as a crystalline solid. A stock solution may be made by dissolving the GW 9662-d₅ in the solvent of choice, which should be purged with an inert gas. GW 9662-d₅ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of GW 9662-d₅ in these solvents is approximately 2, 33, and 35 mg/ml, respectively.

Description

GW 9662 blocks the PPAR γ -induced differentiation of monocytes to osteoclasts by >90% at a dose of 0.1 μ M.¹ It is therefore a much more potent antagonist than BADGE, which is another reported PPAR γ antagonist.¹

Reference

1. Jones, A.E., Wilson, H.K., Meath, P., *et al.* Convenient syntheses of the in vivo carbohydrate metabolites of mycophenolic acid: Reactivity of the acyl glucuronide. *Tetrahedron Lett.* **50(35)**, 4973-4977 (2009).

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, 04/10/2024

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