

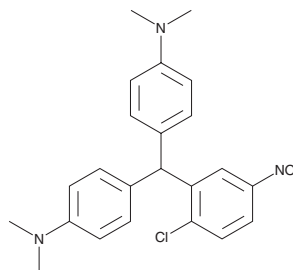
PRODUCT INFORMATION



AL 082D06

Item No. 23455

CAS Registry No.: 256925-03-8
Formal Name: 4,4'-[(2-chloro-5-nitrophenyl)methylene]bis[N,N-dimethyl]-benzenamine
MF: C₂₃H₂₄ClN₃O₂
FW: 409.9
Purity: ≥98%
UV/Vis.: λ_{max}: 268 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

AL 082D06 is supplied as a crystalline solid. A stock solution may be made by dissolving the AL 082D06 in the solvent of choice, which should be purged with an inert gas. AL 082D06 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of AL 082D06 in these solvents is approximately 2.5 and 10 mg/ml, respectively.

AL 082D06 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AL 082D06 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. AL 082D06 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

AL 082D06 is a non-steroidal antagonist of the glucocorticoid receptor (GR; K_i = 210 nM; EC₅₀ = 0.3 nM in a luciferase assay).¹ It is selective for GR over the androgen, progesterone, mineralocorticoid, and estrogen receptors with K_i values greater than 4, 3, 3, and 10 μM, respectively. AL 082D06 inhibits transcriptional activation induced by glucocorticoids and reverses dexamethasone-mediated repression of gene expression in a luciferase reporter assay. It blocks GR binding to DNA *in vitro* when induced by either dexamethasone (Item No. 11015) or mifepristone (Item No. 10006317). In contrast to the steroid-type GR antagonists ZK-299 and mifepristone, GR bound to AL 082D06 does not compete with a constitutively active GR.

Reference

1. Miner, J.N., Tyree, C., Hu, J., *et al.* A nonsteroidal glucocorticoid receptor antagonist. *Mol. Endocrinol.* **17**(1), 117-127 (2003).

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