

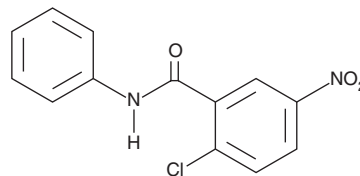
# PRODUCT INFORMATION



## GW 9662

Item No. 70785

**CAS Registry No.:** 22978-25-2  
**Formal Name:** 2-chloro-5-nitrobenzanilide  
**MF:** C<sub>13</sub>H<sub>9</sub>ClN<sub>2</sub>O<sub>3</sub>  
**FW:** 276.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 261 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 is supplied as a crystalline solid. A stock solution may be made by dissolving the GW 9662 in the solvent of choice, which should be purged with an inert gas. GW 9662 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of GW 9662 in these solvents is approximately 2, 33, and 35 mg/ml, respectively.

GW 9662 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GW 9662 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. GW 9662 has a solubility of approximately 0.5 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

GW 9662 blocks the PPAR $\gamma$ -induced differentiation of monocytes to osteoclasts by >90% at a dose of 0.1  $\mu$ M.<sup>1</sup> It is therefore a much more potent antagonist than BADGE, which is another reported PPAR $\gamma$  antagonist.<sup>2</sup>

### References

1. Davies, S.S., Pontsler, A.V., Marathe, G.K., *et al.* Oxidized alkyl phospholipids are specific, high affinity peroxisome proliferator-activated receptor  $\gamma$  ligands and agonists. *J. Biol. Chem.* **276**(19), 16015-16023 (2001).
2. Bendixen, A.C., Shevde, N.K., Dienger, K.M., *et al.* IL-4 inhibits osteoclast formation through a direct action on osteoclast precursors via peroxisome proliferator-activated receptor  $\gamma$ 1. *Proc. Natl. Acad. Sci. USA* **98**, 2443-2448 (2001).

#### 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, 04/08/2024

#### 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