

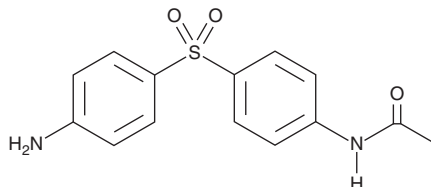
# PRODUCT INFORMATION



## N-acetyl Dapsone

Item No. 17274

**CAS Registry No.:** 565-20-8  
**Formal Name:** N-[4-[(4-aminophenyl)sulfonyl]phenyl]-acetamide  
**Synonym:** NSC 27184  
**MF:** C<sub>14</sub>H<sub>14</sub>N<sub>2</sub>O<sub>3</sub>S  
**FW:** 290.3  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 258, 297, 362 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

N-acetyl Dapsone is supplied as a crystalline solid. A stock solution may be made by dissolving the N-acetyl dapsone in the solvent of choice, which should be purged with an inert gas. N-acetyl Dapsone is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of N-acetyl dapsone in these solvents is approximately 30 mg/ml.

N-acetyl Dapsone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, N-acetyl dapsone should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. N-acetyl Dapsone has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

N-acetyl Dapsone is a metabolite of dapsone, an anti-inflammatory and antibacterial compound that is widely used in the treatment of leprosy, malaria, acne, and various immune disorders.<sup>1</sup> Dapsone is acetylated in the liver to monoacetyldapsone, the major metabolite, and other mono and diacetyl derivatives, and subsequently deacetylated back to diaminodiphenylsulfone (dapsone) until a state of equilibrium is achieved.<sup>1</sup>

### Reference

1. Bonde, S.L., Bhadane, R.P., Gaikwad, A., *et al.* Simultaneous determination of dapsone and its major metabolite N-acetyl dapsone by LC-MS/MS method. *Int. J. Pharm. Pharm. Sci.* **5(Suppl 3)**, 441-446 (2013).

#### 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, 10/14/2022

#### 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