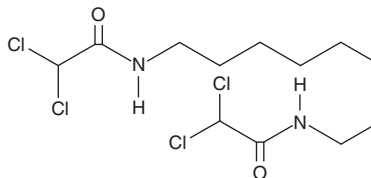


# PRODUCT INFORMATION



**WIN 18,446**  
Item No. 14018

**CAS Registry No.:** 1477-57-2  
**Formal Name:** N,N'-1,8-octanediylbis[2,2-dichloro-acetamide]  
**Synonym:** NSC 59354  
**MF:** C<sub>12</sub>H<sub>20</sub>Cl<sub>4</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 366.1  
**Purity:** ≥98%  
**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

WIN 18,446 is supplied as a crystalline solid. A stock solution may be made by dissolving the WIN 18,446 in the solvent of choice, which should be purged with an inert gas. WIN 18,446 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of WIN 18,446 in these solvents is approximately 0.5, 5, and 10 mg/ml, respectively.

WIN 18,446 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, WIN 18,446 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. WIN 18,446 has a solubility of approximately 0.1 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

Preamble and/or Key Information

## References

1. Reference
2. Reference

**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, 10/13/2022

## 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