

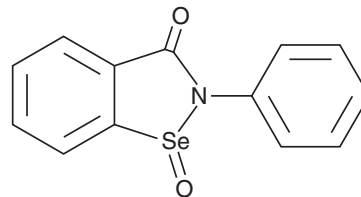
# PRODUCT INFORMATION



## Ebselen Oxide

Item No. 10012298

**CAS Registry No.:** 104473-83-8  
**Formal Name:** 1-oxide-2-phenyl-1,2-benzisoselenazol-3(2H)-one  
**Synonym:** NSC 639772  
**MF:** C<sub>13</sub>H<sub>9</sub>NO<sub>2</sub>Se  
**FW:** 290.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 281 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

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

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of ebselen oxide can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of ebselen oxide in PBS (pH 7.2) is approximately 0.15 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Ebselen is an excellent scavenger of peroxyxynitrite and is used for the treatment of cerebral infarctions. Ebselen oxide, formed by the oxidation of ebselen, lacks antioxidant activity, indicating that it can serve as a negative control for ebselen.<sup>1</sup>

### Reference

1. Lass, A., Witting, P., Stocker, R., *et al.* Inhibition of copper- and peroxy radical-induced LDL lipid oxidation by ebselen: Antioxidant actions in addition to hydroperoxide-reducing activity. *Biochim. Biophys. Acta* **1303(2)**, 111-118 (1996).

#### 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, 04/05/2024

#### 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