

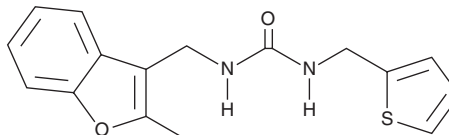
# PRODUCT INFORMATION



**S 07662**

Item No. 45488

**CAS Registry No.:** 883226-64-0  
**Formal Name:** N-[(2-methyl-3-benzofuranyl)methyl]-N'-(2-thienylmethyl)-urea  
**MF:** C<sub>16</sub>H<sub>16</sub>N<sub>2</sub>O<sub>2</sub>S  
**FW:** 300.4  
**Purity:** ≥98%  
**Supplied as:** A 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

S 07662 is supplied as a solid. A stock solution may be made by dissolving the S 07662 in the solvent of choice, which should be purged with an inert gas. S 07662 is soluble (≥10 mg/ml) in DMSO.

## Description

S 07662 is an inverse agonist of the constitutive androstane receptor (CAR).<sup>1</sup> It stabilizes the ligand-binding domain of human CAR, increases the recruitment of nuclear receptor co-repressor (NCOR), and reduces CAR activity (IC<sub>50</sub> = 0.7 μM for the human receptor in a cell-based assay).

## Reference

1. Sengupta, N., Jović, M., Barnaeva, E., et al. A large scale high-throughput screen identifies chemical inhibitors of phosphatidylinositol 4-kinase type II alpha. *J. Lipid Res.* **60(3)**, 683-693 (2019).

### 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, 04/03/2026

## 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