

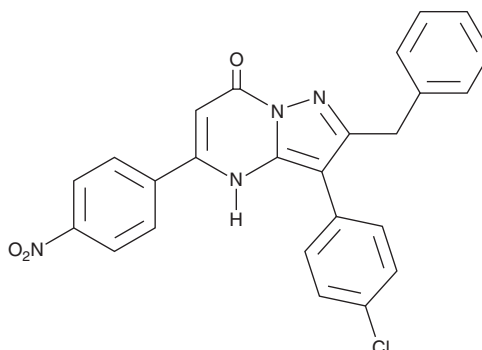
# PRODUCT INFORMATION



## Grassofermata

Item No. 26202

**CAS Registry No.:** 419547-11-8  
**Formal Name:** 3-(4-chlorophenyl)-5-(4-nitrophenyl)-2-(phenylmethyl)-pyrazolo[1,5-a]pyrimidin-7(4H)-one  
**Synonym:** NAV2729  
**MF:** C<sub>25</sub>H<sub>17</sub>ClN<sub>4</sub>O<sub>3</sub>  
**FW:** 456.9  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 272 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

Grassofermata is supplied as a crystalline solid. A stock solution may be made by dissolving the grassofermata in the solvent of choice, which should be purged with an inert gas. Grassofermata is soluble in the organic solvent DMSO at a concentration of approximately 2 mg/ml.

### Description

Grassofermata is an inhibitor of fatty acid transport.<sup>1</sup> It inhibits C<sub>1</sub>-BODIPY-C<sub>12</sub> uptake in HepG2, Caco-2, C2C12, and INS-1E cells (IC<sub>50</sub>s = 8.1-10.6 μM).<sup>1</sup> Grassofermata prevents palmitate-mediated lipid accumulation and cell death in HepG2 cells and primary hepatocytes. *In vivo*, grassofermata (300 mg/kg) decreases absorption of <sup>13</sup>C-oleate in mice.

### Reference

1. Saini, N., Black, P.N., Montefusco, D., *et al.* Fatty acid transport protein-2 inhibitor Grassofermata/CB5 protects cells against lipid accumulation and toxicity. *Biochem. Biophys. Res. Commun.* **465(3)**, 534-541 (2015).

#### 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, 12/21/2022

#### 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