

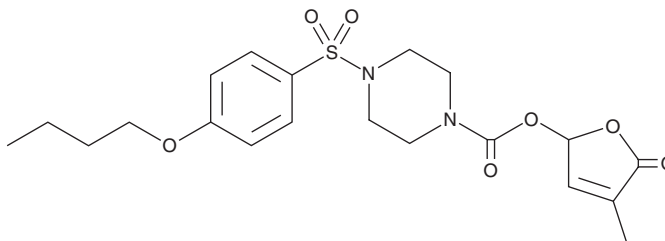
# PRODUCT INFORMATION



## Sphynolactone-7

Item No. 26922

CAS Registry No.: 2305752-57-0  
Formal Name: 4-[(4-butoxyphenyl)sulfonyl]-  
1-piperazinecarboxylic acid,  
2,5-dihydro-4-methyl-5-oxo-2-  
furanylester  
Synonym: SPL7  
MF: C<sub>20</sub>H<sub>26</sub>N<sub>2</sub>O<sub>7</sub>S  
FW: 438.5  
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

Sphynolactone-7 (SPL7) is supplied as a solid. A stock solution may be made by dissolving the SPL7 in the solvent of choice, which should be purged with an inert gas. SPL7 is soluble in DMSO (warmed).

### Description

SPL7 is an agonist of *Striga* hyposensitive to light receptor 7 (ShHTL7), which is found in the parasitic plant *S. hermonthica*.<sup>1</sup> It selectively binds to ShHTL7 (IC<sub>50</sub> = 0.31 μM) over ShHTL2-6, ShHTL9-10, and the strigolactone receptor AtD14 at 10 μM but does inhibit ShHTL8 and ShHTL11 (IC<sub>50</sub>s = 1.2 and 7.8 μM, respectively). SPL7 induces *S. hermonthica* seed germination in the absence of a host, which is fatal to this obligate parasite, with a minimum effective concentration of 10 fM. It induces suicidal germination of *S. hermonthica* seeds and decreases the number of plants emerging from the soil, as well as reduces *S. hermonthica*-induced senescence in maize plants when pre-applied to the soil of co-cultivated maize plants and *S. hermonthica* seeds at a concentration of 10 nM.

### Reference

1. Uraguchi, D., Kuwata, K., Hijikata, Y., *et al.* A femtomolar-range suicide germination stimulant for the parasitic plant *Striga hermonthica*. *Science* **362**(6420), 1301-1305 (2018).

#### 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