

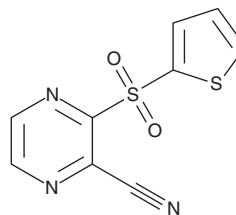
PRODUCT INFORMATION



TSPC

Item No. 35837

CAS Registry No.: 882286-32-0
Formal Name: 3-(2-thienylsulfonyl)-2-pyrazinecarbonitrile
MF: $C_9H_5N_3O_2S_2$
FW: 251.3
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 221, 274 nm
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

TSPC is supplied as a solid. A stock solution may be made by dissolving the TSPC in the solvent of choice, which should be purged with an inert gas. TSPC is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of TSPC in these solvents is approximately 3 mg/ml.

Description

TSPC is an inhibitor of the gibberellin receptor $GID1$.¹ It inhibits the gibberellin-induced interaction between the gibberellin receptor $GID1$ and a negative regulator of its signaling, DELLA, in a yeast two-hybrid assay ($IC_{50} = <30 \mu M$) and gibberellin binding to $GID1$ in the presence or absence of DELLA in a cell-free assay when used at concentrations ranging from 1 to 100 μM . TSPC (10-100 μM) suppresses seed germination and seedling growth in *Arabidopsis*.

Reference

1. Yoon, J.-M., Nakajima, M., Mashiguchi, K., *et al.* Chemical screening of an inhibitor for gibberellin receptors based on a yeast two-hybrid system. *Bioorg. Med. Chem. Lett.* **23**(4), 1096-1098 (2013).

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, 12/15/2022

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