

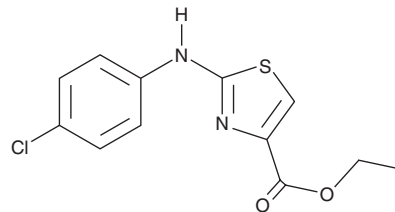
PRODUCT INFORMATION



O412

Item No. 19135

CAS Registry No.: 165682-93-9
Formal Name: 2-[(4-chlorophenyl)amino]-4-thiazolecarboxylic acid, ethyl ester
MF: C₁₂H₁₁ClN₂O₂S
FW: 282.7
Purity: ≥98%
UV/Vis.: λ_{max}: 276 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

O412 is supplied as a crystalline solid. A stock solution may be made by dissolving the O412 in the solvent of choice, which should be purged with an inert gas. O412 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of O412 in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

O412 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, O412 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. O412 has a solubility of approximately 0.12 mg/ml in a 1:7 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Octamer-binding transcription factor 3/4 (Oct3/4) is a transcription factor that has key roles in early embryogenesis and embryonic stem cell pluripotency.^{1,2} O412 is a potent inducer of Oct3/4 expression and translation in human cell lines, including embryonic kidney HEK293, embryonic placental NCCIT, and Oct-deficient HeLa cells.³ It also induces Oct3/4 expression and translation in terminally differentiated human fibroblasts. Through this action, O412 promotes the expression of pluripotency-associated genes in human neonatal foreskin fibroblasts.³

References

1. Niwa, H., Miyazaki, J., and Smith, A.G. Quantitative expression of Oct-3/4 defines differentiation, dedifferentiation or self-renewal of ES cells. *Nat. Genet.* **24**(4), 372-376 (2000).
2. Takahashi, K., Tanabe, K., Ohnuki, M., *et al.* Induction of pluripotent stem cells from adult human fibroblasts by defined factors. *Cell* **131**, 861-872 (2007).
3. Cheng, X., Yoshida, H., Raoofi, D., *et al.* Ethyl 2-((4-Chlorophenyl)amino)thiazole-4-carboxylate and derivatives are potent inducers of Oct3/4. *J. Med. Chem.* **58**(15), 5742-5750 (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, 12/12/2022

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