

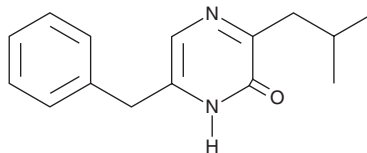
PRODUCT INFORMATION



Pheleuin

Item No. 19269

CAS Registry No.: 169195-23-7
Formal Name: 3-(2-methylpropyl)-6-(phenylmethyl)-2(1H)-pyrazinone
MF: $C_{15}H_{18}N_2O$
FW: 242.3
Purity: $\geq 98\%$
UV/Vis.: λ_{\max} : 324 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

Pheleuin is supplied as a crystalline solid. A stock solution may be made by dissolving the pheleuin in the solvent of choice, which should be purged with an inert gas. Pheleuin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of pheleuin in these solvents is approximately 5 mg/ml in ethanol and 20 mg/ml in DMSO and DMF.

Pheleuin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, pheleuin should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Pheleuin has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Pheleuin is a pyrazinone that has been found in human stool samples.¹ It is produced by gut microbiota nonribosomal peptide synthetases.

Reference

1. Guo, C.-J., Chang, F.-Y., Wyche, T.P., *et al.* Discovery of reactive microbiota-derived metabolites that inhibit host proteases. *Cell* **168**(3), 517-526 (2017).

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, 02/15/2024

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