

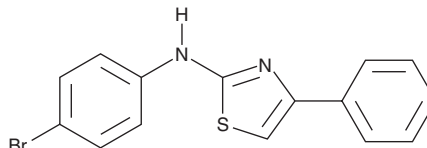
PRODUCT INFORMATION



RCGD 423

Item No. 31420

CAS Registry No.: 108237-91-8
Formal Name: N-(4-bromophenyl)-4-phenyl-2-thiazolamine
MF: $C_{15}H_{11}BrN_2S$
FW: 331.2
Purity: $\geq 98\%$
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

RCGD 423 is supplied as a solid. A stock solution may be made by dissolving the RCGD 423 in the solvent of choice, which should be purged with an inert gas. RCGD 423 is soluble in organic solvents such as ethanol and DMSO. The solubility of RCGD 423 in these solvents is approximately 100 mM.

Description

RCGD 423 is a modulator of gp130.¹ It increases levels of phosphorylated STAT3 and Myc in isolated pig articular chondrocytes ($EC_{50} = 4.5-7.2 \mu M$), effects that can be reversed by the gp130 inhibitor SC144 (Item No. 29714). RCGD 423 reduces cellular apoptosis and hypertrophy induced by bone morphogenic protein 4 (BMP4) in isolated mouse total limb cells and isolated human healthy and osteoarthritic articular chondrocytes. *In vivo*, RCGD 423 (4 μg /joint) reduces cartilage loss in a rat model of osteochondral injury.

Reference

- Shkyhan, R., Handel, B.V., Bogdanov, J., *et al.* Drug-induced modulation of gp130 signalling prevents articular cartilage degeneration and promotes repair. *Ann. Rheum. Dis.* **77**(5), 760-769 (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, 11/09/2022

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