

PRODUCT INFORMATION



Kobe 2602

Item No. 16262

CAS Registry No.: 454453-49-7

Formal Name: 2-[2,6-dinitro-4-(trifluoromethyl)phenyl]-N-(4-fluorophenyl)-hydrazinecarbothioamide

MF: $C_{14}H_9F_4N_5O_4S$

FW: 419.3

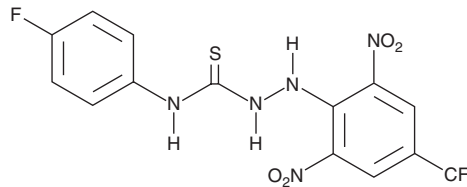
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 226, 232, 366 nm

Supplied as: A crystalline 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

Kobe 2602 is supplied as a crystalline solid. A stock solution may be made by dissolving the kobe 2602 in the solvent of choice, which should be purged with an inert gas. Kobe 2602 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of kobe 2602 in these solvents is approximately 5, 20, and 10 mg/ml, respectively.

Kobe 2602 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, kobe 2602 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Kobe 2602 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

The Ras family of small GTPases (H-Ras, K-Ras, and N-Ras) function as molecular switches, cycling between a GTP-bound active state and a GDP-bound inactive state, to turn on downstream Raf protein kinases. This initiates complex signaling pathways involved in cell growth, differentiation, and apoptosis. Mutations leading to aberrant Ras activation are frequently associated with various human cancers. Kobe 2602 is a selective Ras inhibitor that blocks H-Ras GTP binding to c-Raf-1 ($K_i = 149 \mu M$).¹ Kobe 2602 has been shown to inhibit both anchorage-dependent and -independent growth and to induce apoptosis of H-Ras^{G12V}-transformed NIH 3T3 cells ($IC_{50} = 1.4-2 \mu M$).¹ At an oral dose of 80 mg/kg, it also exhibits antitumor activity in mice bearing a xenograft of human colon cancer SW480 cells expressing K-Ras^{G12V}.¹

Reference

1. Shima, F., Yoshikawak, Y., Ye, M., *et al.* In silico discovery of small-molecule Ras inhibitors that display antitumor activity by blocking the Ras-effector interaction. *Proc. Natl. Acad. Sci. USA* **110**(20), 8182-8187 (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, 10/28/2022

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