

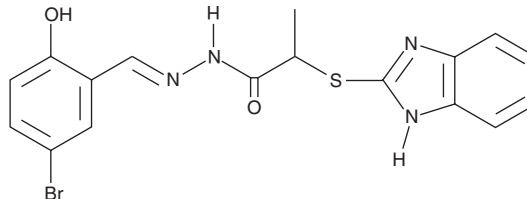
PRODUCT INFORMATION



KH7

Item No. 13243

CAS Registry No.: 330676-02-3
Formal Name: 2-(1H-benzimidazol-2-ylthio)-2-[(5-bromo-2-hydroxyphenyl)methylene]hydrazide, propanoic acid
MF: C₁₇H₁₅BrN₄O₂S
FW: 419.3
Purity: ≥98%
UV/Vis.: λ_{max}: 284, 337 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

KH7 is supplied as a crystalline solid. A stock solution may be made by dissolving the KH7 in the solvent of choice, which should be purged with an inert gas. KH7 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of KH7 in these solvents is approximately 0.5, 25, and 30 mg/ml, respectively.

Description

Soluble adenylyl cyclase mediates bicarbonate-stimulated production of the ubiquitous second messenger adenosine 3',5'-cyclic mononucleotide (cAMP).¹ It is abundantly expressed in sperm, in fluid transporting tissues such as kidney cortex and medulla, and in other bicarbonate-responsive tissues and cells. KH7 is a selective inhibitor of soluble adenylyl cyclase (sAC) that has little effect on transmembrane adenylyl cyclases.² It displays an IC₅₀ value between 3-10 μM toward sAC.² At 50 μM, KH7 decreases basal cAMP accumulation in sperm.² KH7 has been used in diverse research applications including cAMP-mediated signaling events required for mammalian egg fertilization, sAC regulation of Na⁺ transport in the kidney, and mitochondrial-dependent apoptosis in response to various stress stimuli.²⁻⁴ KH7 is not active against sAC in the presence of detergents. At concentrations above 50 μM, KH7 exhibits non-specific membrane disruption effects when used on cells.

References

1. Wuttke, M.S., Buck, J., and Levin, L.R. Bicarbonate-regulated soluble adenylyl cyclase. *JOP* **2(4 Suppl)**, 154-158 (2001).
2. Hess, K.C., Jones, B.H., Marquez, B., et al. The "soluble" adenylyl cyclase in sperm mediates multiple signaling events required for fertilization. *Dev. Cell* **9(2)**, 249-259 (2005).
3. Hallows, K.R., Wang, H., Edinger, R.S., et al. Regulation of epithelial Na⁺ transport by soluble adenylyl cyclase in kidney collecting duct cells. *J. Biol. Chem.* **284(9)**, 5774-5783 (2009).
4. Kumar, S., Kostin, S., Flacke, J.P., et al. Soluble adenylyl cyclase controls mitochondria-dependent apoptosis in coronary endothelial cells. *J. Biol. Chem.* **284(22)**, 14760-14768 (2009).

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, 03/11/2024

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