

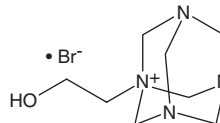
PRODUCT INFORMATION



Y11

Item No. 18521

CAS Registry No.: 1086639-59-9
Formal Name: 1-(2-hydroxyethyl)-3,5,7-triaza-1-azoniatricyclo[3.3.1.1^{3,7}]decane, monobromide
Synonyms: FAK Inhibitor 10, NSC 206142
MF: C₈H₁₇BrN₄O
FW: 265.2
Purity: ≥95%
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

Y11 is supplied as a crystalline solid. A stock solution may be made by dissolving the Y11 in the solvent of choice, which should be purged with an inert gas. Y11 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of Y11 in these solvents is approximately 10 and 0.3 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of Y11 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of Y11 in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Focal adhesion kinases (FAKs) are non-receptor tyrosine kinases that play roles in regulating diverse processes, including cell adhesion, spreading, migration, proliferation, and apoptosis. They are over-expressed in many types of cancer. Y11 inhibits FAK1 autophosphorylation by blocking phosphorylation of Y397.¹ At 100 μM, it was shown to decrease cell viability of T47D breast cancer and C8161 melanoma cell lines.¹

Reference

1. Golubovskaya, V.M., Nyberg, C., Zheng, M., *et al.* A small molecule inhibitor, 1,2,4,5-benzenetetraamine tetrahydrochloride, targeting the Y397 site of focal adhesion kinase decreases tumor growth. *J. Med. Chem.* **51**(23), 7405-7416 (2008).

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, 11/01/2022

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