

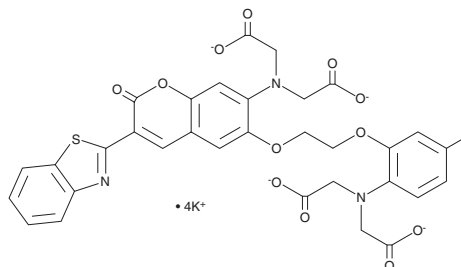
PRODUCT INFORMATION



BTC (potassium salt)

Item No. 20423

CAS Registry No.: 216453-54-2
Formal Name: N-[3-(2-benzothiazolyl)-6-[2-[2-bis(carboxymethyl)amino]-5-methylphenoxy]ethoxy]-2-oxo-2H-1-benzopyran-7-yl]-N-(carboxymethyl)glycine, tetrapotassium salt
MF: C₃₃H₂₅N₃O₁₂S • 4K
FW: 844.0
Purity: ≥90%
Supplied as: A 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

BTC (potassium salt) is supplied as a solid. A stock solution may be made by dissolving the BTC (potassium salt) in the solvent of choice. BTC (potassium salt) is soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

BTC is a low affinity calcium indicator ($K_d = 7-26 \mu\text{M}$) that displays excitation/emission spectra of 401/529 nm, respectively. It exhibits a shift in the excitation maximum from approximately 480 to 401 nm upon calcium binding, enabling determination of ratiometric calcium measurements.¹ BTC is suitable for detecting elevated calcium levels associated with activation of smooth muscle, neurons, and intracellular calcium stores.^{1,2}

References

1. Sabnis, R.W. Handbook of biological dyes and stains: Synthesis and industrial applications. John Wiley & Sons, Inc., Hoboken, NJ, USA (2010).
2. Hyrc, K.L., Bownik, J.M., and Goldberg, M.P. Neuronal free calcium measurement using BTC/AM, a low affinity calcium indicator. *Cell Calcium* 24(3), 165-175 (1998).

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, 12/19/2022

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