

PRODUCT INFORMATION



IPG-4 TMA

Item No. 35539

Synonyms: APG-4, Asante Potassium Green-4,
Ion Potassium Green-4

MF: $C_{58}H_{64}Cl_2N_3O_{15} \cdot 3C_4H_{12}N$

FW: 1336.5

Purity: $\geq 85\%$

UV/Vis.: λ_{max} : 524 ± 3 nm

Ex./Em. Max: 525/545 nm

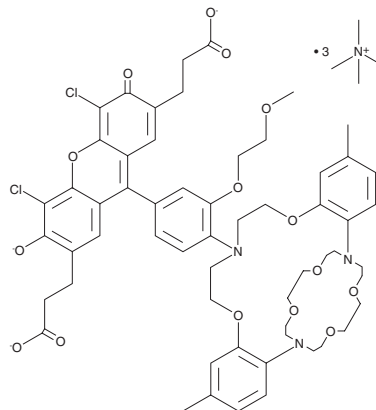
Supplied as: A solid

Storage: 20°C

Stability: ≥ 4 years

Special Conditions: Protect from light and moisture

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.



Description

IPG-4 TMA is a cell-impermeable fluorescent potassium indicator that has been used to detect extracellular potassium in *B. subtilis* biofilms.¹ It binds to potassium ($K_d = 7$ mM) and displays excitation/emission maxima of 525/545 nm, respectively. IPG-4 TMA is also available in a cell-permeable form (Item No. 35541).

Reference

1. Prindle, A., Liu, J., Asally, M., *et al.* Ion channels enable electrical communication in bacterial communities. *Nature* **527(7576)**, 59-63 (2015).

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, 12/02/2022

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