

PRODUCT INFORMATION



IPG-4 AM

Item No. 35541

CAS Registry No.: 3019514-72-5

Synonyms: APG-4 Acetoxymethyl ester,
APG-4 AM Asante Potassium Green-4
Acetoxymethyl ester, Asante Potassium Green-4 AM,
Ion Potassium Green-4 Acetoxymethyl ester,
IPG-4 Acetoxymethyl ester

MF: $C_{67}H_{79}Cl_2N_3O_{21}$

FW: 1,333.3

Purity: $\geq 90\%$

UV/Vis.: λ_{max} : 524 nm

Ex./Em. Max: 525/545 nm

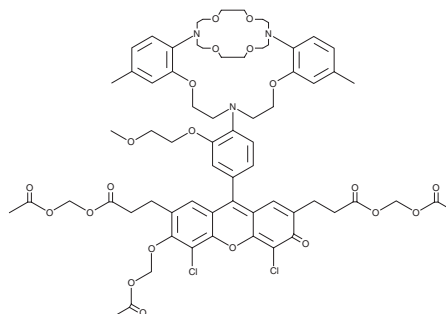
Supplied as: A solid

Storage: $-20^{\circ}C$

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.



Laboratory Procedures

IPG-4 AM is supplied as a solid. A stock solution may be made by dissolving the IPG-4 AM in the solvent of choice, which should be purged with an inert gas. IPG-4 AM is soluble in DMSO.

Description

IPG-4 AM is a membrane-permeable fluorescent potassium indicator that has been used to detect intracellular potassium.¹ It binds to potassium ($K_d = 7$ mM) and displays excitation/emission maxima of 525/545 nm, respectively. IPG-4 AM is also available in a cell-impermeable form (Item No. 35539).

Reference

1. Rana, P.S., Gibbons, B.A., Vereninov, A.A., *et al.* Calibration and characterization of intracellular Asante Potassium Green probes, APG-2 and APG-4. *Anal. Biochem.* **567**, 8-13 (2019).

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, 03/29/2024

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