

PRODUCT INFORMATION



IPG-2 AM

Item No. 35540

CAS Registry No.: 1369302-24-8
Formal Name: 6-[(acetyloxy)methoxy]-4,5-dichloro-9-[3-methoxy-4-[16-(2-methoxy-4-methylphenyl)-1,4,10,13-tetraoxa-7,16-diazacyclooctadec-7-yl]phenyl]-3-oxo-3H-xanthene-2,7-dipropionic acid, 2,7-bis[(acetyloxy)methyl] ester

Synonyms: APG-2 Acetoxymethyl ester, APG-2 AM, Asante Potassium Green-2 Acetoxymethyl ester, Asante Potassium Green-2 AM, ION Potassium Green-2 Acetoxymethyl ester, ION Potassium Green-2 AM, IPG-2 Acetoxymethyl ester

MF: $C_{55}H_{64}Cl_2N_2O_{19}$

FW: 1,128.0

Purity: $\geq 90\%$

Ex./Em. Max: 525/545 nm

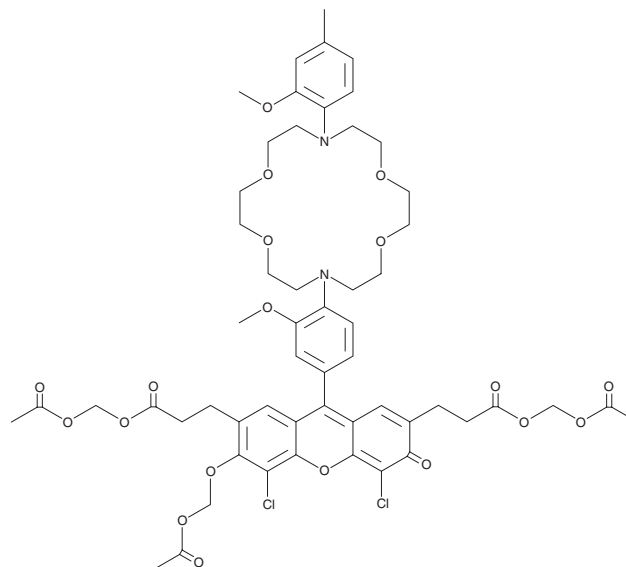
Supplied as: A solid

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.



Description

IPG-2 AM is a cell-permeable fluorescent potassium indicator. It binds to potassium ($K_d = 18$ mM) and displays excitation/emission maxima of 525/545 nm, respectively. IPG-2 AM has been used to kinetically monitor potassium levels in isolated human platelets via flow cytometry.¹ IPG-2 is also available in a cell-impermeable form (Item No. 35533).

Assay Protocol

Note: Allow all reagents to warm to room temperature before proceeding.

1. Add 10 ml of assay buffer to a 15 or 50 ml conical tube.

Note 1: HEPES-buffered Hank's balanced salt solution (HBSS), pH 7.2-7.4, is recommended, although other buffers can be used.

2. Add 100 μ l of a 100X Pluronic™ F-127 solution (1-50% w/v) to the conical tube*. Pluronic™ F-127 is a biocompatible surfactant used to ensure equitable dye distribution and cellular loading.
 - a. Optional: Add 100 μ l of 2 mM probenecid stock solution to the conical tube. Probenecid (Item No. 14981) is an anion transport inhibitor used to improve intracellular dye retention. Use of probenecid is recommended, but not required, for all cell types and dyes.

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, 12/12/2022

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

PRODUCT INFORMATION



**Final working concentration of Pluronic™ F-127 should be between 0.01 and 0.5% w/v. User should optimize the concentration of Pluronic™ F-127 to suit experimental requirements.*

Note 2: Probenecid is an inhibitor or agonist of multiple ion channels and may have undesirable cellular effects that could affect dye performance.

3. Vortex conical tube briefly to mix.
4. Dissolve IPG-2 AM in 25 μ l of DMSO and vortex dye tube briefly to mix.
5. Centrifuge dye tube briefly to collect all contents at the tube bottom.
6. Add entire contents of dye tube to the conical tube containing the assay buffer solution to make the dye loading solution.
7. Vortex conical tube briefly to mix.

Note 3: The dye loading solution should be used within two hours for best results.

8. Remove cell culture medium and add dye loading solution. Recommended volumes are:
 - a. 35 mm dish or 6-well plate: 1.5 ml/dish or well
 - b. 96-well plate: 100 μ l/well
 - c. 384-well plate: 20 μ l/well

Note 4: To prevent cell detachment or if using suspension cells, the dye loading solution can be added directly to the media-containing wells. User must double the component concentrations to achieve the same final concentration of all reagents.

9. Incubate cells with the dye loading solution at 37°C for 60 minutes.
10. Read fluorescence using a plate reader at excitation and emission wavelengths of 525 and 545 nm, respectively.
Or
Image using a fluorescence microscope with filters for YFP, GFP, or fluorescein.

Reference

1. Aliotta, A., Calderara, D.B., and Alberio, L. Flow cytometric monitoring of dynamic cytosolic calcium, sodium, and potassium fluxes following platelet activation. *Cytometry A*. **97(9)**, 933-944 (2020).

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