

# Rhod-2 AM (special packaging)

PromoKine

-[2-Amino-5-(3-dimethylamino-6-dimethylammonio-9-xanthenyl)phenoxy]-2-(2-amino-5-methylphenoxy)ethane-N,N,N',N'-tetraacetic acid, tetraacetoxymethyl ester

## Instruction Manual

Catalog Number	PK-CA707-50023
Description	<p>Rhod-2 has the longest fluorescent wavelength signal of all the calcium indicators. It has a rhodamine-like fluorophore whose excitation and emission maxima are 556 nm and 576 nm, respectively. This makes it a convenient excitation source for argon and krypton lasers. Although it was thought that the fluorescent signal of Rhod-2 only increases several times with the calcium complex, Rhod-2 increases its signal with calcium 80-100 times because of its high purity. Its signal intensity is the strongest of all the calcium probes. Rhod-2 is thus highly recommended as a probe for intracellular calcium monitoring using laser microscopes. Rhod-2 is reported to have a better loading profile at the point of localization, especially in neural slice cultures. The dissociation constant of Rhod-2 with calcium (<math>K_d=1.0</math> mM) is the highest of all the fluorescent calcium probes, providing a wider range for monitoring calcium concentration. Rhod-2 is similar to Fluo-3 in that the excitation and emission spectra do not undergo a shift and the sensor is essentially nonfluorescent before <math>Ca^{2+}</math> binding but becomes more fluorescent with increasing <math>Ca^{2+}</math> concentration. The absorption and emission maxima of Rhod-2, however, are longer than those of Fluo-3. The longer absorption and emission wavelengths of Rhod-2 may make it useful for some applications where autofluorescence is a problem, or where another fluorescent dye of shorter wavelengths is used at the same time. The fluorescent enhancement for Rhod-2 from low <math>[Ca^{2+}]</math> to high <math>[Ca^{2+}]</math> was smaller than that for Fluo-3, and also in general Rhod-2 is somewhat less fluorescent than Fluo-3.</p> <p>Rhod-2 AM is an acetoxymethyl ester derivative of Rhod-2 that can be easily loaded into cells via incubation using the AM method. Because of the relatively low water solubility of the AM ester, Pluronic F-127 (See PK-CA707-59000 and refs. therein), a mild detergent, is often used as a dispersing agent to facilitate the loading. Rhod-2/AM itself does not bind <math>Ca^{2+}</math>, but it is readily hydrolyzed to Rhod-2 by endogenous esterases once the dye is inside the cells. For other information on Rhod-2, see PK-CA707-50020/50021/50022.</p> <p>PromoKine offers A-23187 (PK-CA707-59001), an ionophore that is commonly used for intracellular calibration of calcium indicators. PromoKine also offers EDC (PK-CA707-59002, also known as EDAC), which can be used to fix calcium indicators in cells, if post histochemical studies are desired following physiological experiments.</p>
Quantity	10 x 100 $\mu$ g
Excitation / Emission Maxima	$\lambda_{exc} \backslash \lambda_{em}$ (DNA) = 556/576 nm (low or high $[Ca^{2+}]$ ); Extinction Coefficient: 80,000 $M^{-1}cm^{-1}$ (556 nm)
Molecular Structure	
Molecular Weight / Molecular Formula	1132 Da; $C_{52}H_{59}BrN_4O_{19}$
Purity	>95% (as determined by TLC)
Appearance / Formulation / Solubility	Red solid; soluble in DMSO
Intended Use	For in vitro research use only. Not for diagnostic or therapeutic procedures.
Storage & Stability	Stored desiccated at $-20^{\circ}C$ . Protect from light, especially when in solution.

<b>Applications</b>	Fluorescent calcium indicator; see Description
<b>References</b>	1) Meyer, T., et al. Biochemistry 29, 32(1990) 2) Minta, A., et al. J. Biol. Chem. 264, 8171(1989).
<b>Caution</b>	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

**PromoCell GmbH**

Sickingenstr. 63/65  
69126 Heidelberg  
Germany

Email: [info@promokine.info](mailto:info@promokine.info)  
[www.promokine.info](http://www.promokine.info)

**North America**

Phone: 1 - 866 - 251 - 2860 (toll free)  
Fax: 1 - 866 - 827 - 9219 (toll free)

**Deutschland**

Telefon: 0800 - 776 66 23 (gebührenfrei)  
Fax: 0800 - 100 83 06 (gebührenfrei)

**France**

Téléphone: 0800 90 93 32 (ligne verte)  
Téléfax: 0800 90 27 36 (ligne verte)

**United Kingdom**

Phone: 0800 - 96 03 33 (toll free)  
Fax: 0800 - 169 85 54 (toll free)

**Other Countries**

Phone: +49 6221 - 649 34 0  
Fax: +49 6221 - 649 34 40