

Mag-Fura-2 tetrapotassium salt *Cell impermeant*

Catalog number: 20380

Unit size: 1 mg

Product Details

Storage Conditions Freeze (<-15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

Chemical Properties

Appearance Yellow solid

Molecular Weight 586.68

Soluble In Water

Chemical Structure

Spectral Properties

Excitation Wavelength 336 nm

Emission Wavelength 505 nm

Applications

Mag-Fura-2, tetrapotassium salt is an intracellular magnesium indicator that is ratiometric and UV light-excitable. This water-soluble salt form is useful for intracellular loading by microinjection or infusion from patch pipette. This particular magnesium indicator is also used for high concentration calcium detection.