

Product Information Sheet

Ordering Information

Product Number:	20404
Product Name:	Mag-520™ potassium salt
Unit Size:	10x50 ug
Storage Conditions:	Freeze (<-15 °C), Minimize light exposure
Expiration Date:	12 months upon receiving

Chemical and Spectral Properties

Appearance:	Solid
Molecular Weight:	475.45
Soluble In:	Water
Excitation Wavelength:	N/A
Emission Wavelength:	N/A

Application Notes

A vast majority of the existing magnesium ion indicators are based on tricarboxylate APTRA chelator derived from the popular tetracarboxylate BAPTA calcium ion chelator. They include mag-fura-2, mag-indo-1, mag-fluo-4 and mag-rhod-3. However, all of them have higher affinity for calcium than magnesium although they were designed for detecting magnesium ion. Mag-520[™] is the first commercial magnesium indicator that has higher affinity for magnesium than calcium. Its significantly improved selectivity can be used for measuring magnesium ion with minimal interference from calcium ion compared to other commercial magnesium ion indicators such as the popular mag-fluo-4.