

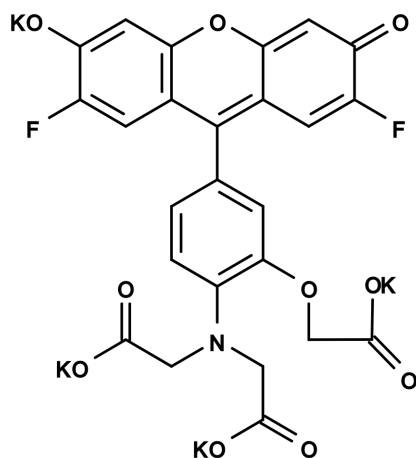
Product Information Sheet

Ordering Information

| | |
|---------------------|---|
| Product Number: | 20400 |
| Product Name: | Mag-Fluo-4 potassium salt |
| Unit Size: | 10x50 ug |
| Storage Conditions: | Freeze (<-15 °C), Desiccated, Avoid Light |
| Expiration Date: | 12 months upon receiving |

Chemical and Spectral Properties

| | |
|---------------------|--------|
| Appearance: | Solid |
| Molecular Weight: | 681.77 |
| Chemical Structure: | |



| | |
|------------------------|-------|
| Soluble In: | Water |
| Excitation Wavelength: | 494 |
| Emission Wavelength: | 516 |

Application Notes

The cell-impermeant Mag-Fluo-4 is an analog of Fluo-4 with a K_d for Mg ion of 4.7 mM and a K_d for Ca ion of 22 μ M, making it useful as an intracellular Mg ion indicator as well as a low-affinity Ca ion indicator.