

Product Information Sheet

Ordering Information

Product Number: 1375
Product Name: iFluor™ 810 acid
Unit Size: 1 mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 24 months upon receiving

Chemical, Physical and Spectral Properties

Molecular Weight: 1373.43
Appearance: Blue/Green powder
Soluble in: Water, DMF or DMSO
Excitation Wavelength: 809 nm
Emission Wavelength: 821 nm

Application Notes

In vivo fluorescence imaging uses a sensitive camera to detect fluorescence emission from fluorophores in whole-body living small animals. To overcome the photon attenuation in living tissue, fluorophores with long emission at the infrared (IR) region are generally preferred. Recent advances in imaging strategies and reporter techniques for in vivo fluorescence imaging include novel approaches to improve the specificity and affinity of the probes and to modulate and amplify the signal at target sites for enhanced sensitivity. Further emerging developments are aiming to achieve high-resolution, multimodality and lifetime-based in vivo fluorescence imaging. Our iFluor™ 810 is designed to label proteins and other biomolecules with infrared fluorescence. Conjugates prepared with iFluor™ 810 have the excitation and emission in the IR range. iFluor™ 810 dye emission is well separated from commonly used far-red fluorophores such as Cy5, Cy7 or allophycocyanin (APC), facilitating multicolor analysis. This fluorophore is also useful for small animal in-vivo imaging applications or for other imaging applications that require IR detections.