

# Congo Red NHS Ester

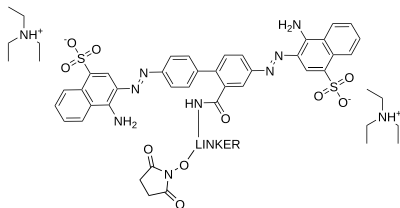
Catalog Number: 23069

Unit Size: 5 mg

## Product Details

|                    |  |
|--------------------|--|
| Storage Conditions | Freeze (< -15 °C), Minimize light exposure |
| Expiration Date    | 12 months upon receiving                   |

## Chemical Properties

|                    |  |
|--------------------|--|
| Appearance         | Solid red  |
| Molecular Weight   | 1109.33  |
| Soluble In         | DMSO   |
| Chemical Structure |  |

## Spectral Properties

|                       |     |
|-----------------------|-----|
| Excitation Wavelength | N/A |
| Emission Wavelength   | N/A |

## Applications

Congo Red has been used as a common qualitative staining method for the identification of amyloids in vitro and in tissue sections. Congo Red NHS ester is an excellent building block to develop Congo Red-based amyloid probes. The investigation of the amyloid formation is currently one of the most important tasks in the neurological field. Research in this area involves not only understanding how amyloidosis develops and how to prevent amyloid plaque formation but also investigating functional amyloids. In amyloidosis diseases, amyloids can be detected in the kidney, liver, brain and other human and animal organs. The development of neurodegenerative human diseases, such as Alzheimer's disease and Parkinson's disease, is probably the consequence of accumulation of insoluble amyloid plaques in the nerve tissue.