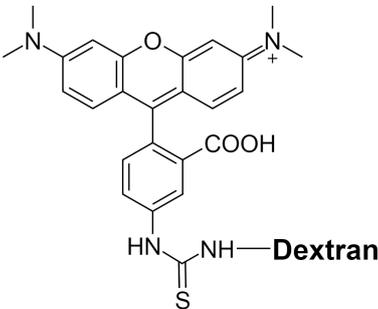
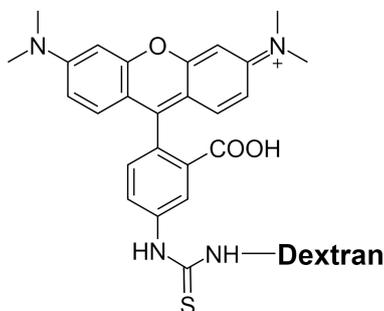


TRITC–dextran conjugate (average MW = ~150K)Catalog number: 21710
Unit size: 25 mg**Product Details**

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

Chemical Properties

Appearance	Solid
Molecular Weight	N/A
Soluble In	Water
Chemical Structure	

**Spectral Properties**

Excitation Wavelength	544 nm
Emission Wavelength	570 nm

Applications

TRITC-labeled dextran is used in cardiovascular, microcirculation, perfusion, cell monolayer and cell membrane permeability research as fluorescent flux tracer compound that supports the measurement of processes such as blood flow, membrane damage, vascular drainage and renal elimination. It is also used in microcirculation and cell permeability research utilizing microfluorimetry, and possible in perfusion studies in animals. TRITC-dextran has also been used to study plant cell wall porosity and capillary permeability. Plasma proteins have been shown not to bind to TRITC-dextran. TRITC-dextran of 150 kDa may be used to help assess the integrity of the blood brain barrier (BBB) and to perform fluorescence microlymphography. It may also be used as a microcirculation vascular flux tracer.