



Lectin Kit I, Fluorescein FLK-2100

Product Images



Short Description

Our lectin screening kits are designed to provide the investigator with a panel of seven lectins or lectin conjugates. The lectins have been selected to offer a variety of sugar specificities and are of the same high quality as the reagents offered individually.

Each Kit I (BK-1000, FLK-2100, RLK-2200) consists of 1 mg of the following lectins or lectin conjugates: Con A, DBA, PNA, RCA120, SBA, UEA I, WGA.

Additional Information

Unit Size	1 kit
Applications	Immunofluorescence, Glycobiology
Conjugate	Fluorescein
Color of Fluorescence	Green
Sugar Specificity	Mannose, Galactose, Fucose, N-Acetylgalactosamine, N-Acetylglucosamine

