



Lectin Kit II, Biotinylated

Product Images



Short Description

Our lectin screening kits are designed to provide the investigator with a panel of six 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 II (LK-3000, BK-2100, FLK-3100, RLK-3200) consists of 1 mg of the following lectins or lectin conjugates: GSL I, LCA, PHA-E, PHA-L, PSA, Succinylated WGA.

Additional Information

Unit Size	1 kit
Applications	Immunohistochemistry / Immunocytochemistry, Immunofluorescence, Blotting Applications, Elispot, ELISAs, Glycobiology
Conjugate	Biotinylated

