



CERTIFICATE OF ANALYSIS

Product FLUORESCEIN PEANUT AGGLUTININ

Catalog No. EL-1071

Amount 10 mg

Lot No. ZC1102

Concentration 5 mg/ml

Solution 10 mM HEPES, 0.15 M NaCl, pH 7.5, 0.08% sodium azide,
0.1 mM Ca⁺⁺

Free Fluorescein None

F/P (molar) 4.2

Maximum Excitation @ 495-500 nm

Maximum Emission @ 514-521 nm

Active Conjugate 99%

Storage Conditions Refrigerate in the dark

Other Remarks:

A precipitate may form during storage. This does not have a significant adverse effect on the product. If a precipitate forms upon long-term storage at 4 °C, warm to 37 °C and centrifuge before use.

The recommended concentration range for use is 5-20 µg/ml.

Inhibiting/Eluting Sugar: 200 mM galactose (S-9003).

For immunofluorescence applications, we recommend that sections be mounted in VECTASHIELD® Hard♦Set™ Mounting Medium (H-1400) or VECTASHIELD® Hard♦Set™ Mounting Medium with DAPI (H-1500).