



CERTIFICATE OF ANALYSIS

Product **FLUORESCEIN PEANUT AGGLUTININ**

Catalog No. **EL-1071**

Amount **5 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⁺⁺, 0.01 mM Mn⁺⁺**

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).