



CERTIFICATE OF ANALYSIS

Product FLUORESCHEIN SAMBUCUS NIGRA (ELDERBERRY) BARK LECTIN

Catalog No. EL-1301

Amount 2 mg

Lot No. ZB0106

Concentration 2 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) 7.7

Active Conjugate 95%

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.