



CERTIFICATE OF ANALYSIS

Peridinin-Chlorophyll Protein Complex (PerCP)

Product Number D11-000
Lot Number XXXX
Approximately 20mg/ml
Concentration 23.1 mg/mL
Store at 2-8°C

Form/Shipping

Supplied as a liquid, store at 2-8° C in the dark. Do not freeze. PerCP is sensitive to freeze-thaw cycle.

Handling

Avoid exposure to light and heat. Please use an AU/mg of 11.6 to determine PerCP concentration.

Buffer

50 mM HEPES (pH 7.5) and 15 mM sodium azide

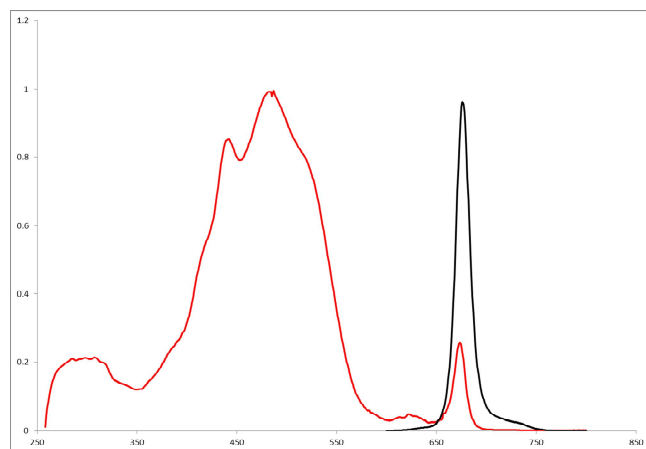
Note

For research use only, not for diagnostic or therapeutic use

<u>Specifications</u>	<u>Actual</u>
Absorbance Max (483 +/- 1nm)	482 nm
Emission Max (676 +/-1 nm)	676.5 nm
A ₄₈₃ /A ₂₈₀ (4.25)	4.78
HPLC-SEC	100.0%

Method

The PerCP concentration was measured in 50 mM HEPES (pH 7.5) + 15 mM sodium azide. The reported results are the average of three different measurements.



Absorbance and emission spectra of Peridinin- Chlorophyll Protein Complex (PerCP) in 10 mM Sodium Phosphate (pH 7.4). Emission scan was taken with an excitation at 488nm.