
RayBiotech Searchlight Ultra-Sensitive Chemiluminescent HRP Membrane/Blotting Substrate

(Two Components A + B)

Components

The substrate kit is supplied as two components: Reagent A (Part #. HRP-A), Reagent B (Part #. HRP-B).

- Catalog #. RB-SRCHLGT-200 (200 ml: 100 ml Reagent A, 100 ml Reagent B)
- Catalog #. RB-SRCHLGT-50 (50 ml: 25 ml Reagent A, 25 ml Reagent B)

Product Description

RayBiotech Searchlight Ultra-Sensitive Chemiluminescent HRP Membrane/Blotting Substrate kit can be used for high-sensitivity detection of horseradish peroxidase (HRP) in membrane/dot blotting and Western blotting. Raybiotech chemiluminescent substrates provide superior sensitivity compared to competitor products and are compatible with X-ray film and digital CCD imaging.

Storage

2-8 °C. Protected from light and heat. Keep container tightly closed.

Stability

Stable for 24 months when stored at 2-8 °C and protected from light.

Use Instruction

The chemiluminescent substrate should be prepared immediately before use. In a clean container, mix one part of Reagent A with one part of Reagent B to a total volume appropriate to your application (we recommend: use 50-100 µl mixed substrate per cm² membrane). Mix well and protect the substrate from light while using. The diluted mixture of Reagents A and B is stable for several hours in the dark.

MSDS & COA

Please download from www.raybiotech.com.

**The products are furnished for LABORATORY RESEARCH USE ONLY.
Not for therapeutic use.**