

HTRF® compatibility certificate

Upon the results of the assessment procedure, Cisbio/Perkin Elmer certifies that Berthold's LB 942 TriStar 5 plate reader is approved for HTRF assay readout.

- Certifying tests

Europium donor / Red acceptor combination	✓ Validated
Terbium donor / Red acceptor combination	✓ Validated

- Color plate restriction

White and Black plates	White plates only
------------------------	-------------------

- Optional tests (not used for the certification)

Terbium donor / Green acceptor combination	✓ Validated
--	-------------

Codolet, August 11th, 2022

Sandrine Cabirol
Instrument Certification Officier



Delphine Jaga
Research Reagent Application SME Leader
Instrument Certification Supervisor


