

MANUAL

Technology: HTRF™

Protein-Protein Interaction

d2 labeling kit

Storage details		Storage upon receipt
Set sent in dry ice	8 nmoles d2 dye	-60°C or below
Set sent at room temperature (RT)	Buffers and column	2-8°C

Part number: 62D2DPEA

Version: 06

Date: July 2024

KIT DESCRIPTION AND INTENDED USE

The HTRF d2 labeling kit is a comprehensive set of reagents that enables the covalent coupling of d2 acceptor dye to biomolecules.

The kit is suitable for labeling antibodies and proteins of similar molecular weights, N-hydroxysuccinimide-activated d2 dye reacts with primary amines to form stable dye-conjugates under mild conditions.

d2 acceptor, an organic motif of approximately 800 Da, is highly compatible with HTRF cryptates. Its excitation spectrum overlaps that of Eu₃⁺ and Lumi4™-Tb cryptate emission, allowing the donor to excite d2 with a high quantum yield. d2's maximum light output is at 665nm.

Note: The kit was designed as a tool for assessing the biological activity of d2-labeled moieties. For any enquiry about custom labeling services, please directly contact Revvity.

EQUIPMENT AND MATERIALS REQUIRED BUT NOT INCLUDED

- Precision micropipettes with disposable tips, capable of dispensing 10-1000 µL.
- Column stand and clamp.
- Vortex.
- Test tubes.
- HTRF™ compatible reader (more information about compatible reader on our website)

SUPPLIED REAGENTS AND STABILITY

Supplied reagents

d2 labeling reagent 8 nmoles (6µg) - desiccated	1 microtube + desiccant	Store at -60°C or below until use
Purification column	1 column	Store at 2-8°C
Elution buffer (100 mM phosphate buffer pH7, 2 mM NaN ₃)	1 vial, 20 mL	Store at 2-8°C until use

Reagent stability

The d2 dye should be stored at -60°C or below, the other 2 components must be stored at 2-8°C before use. Under these conditions, the kit is stable until the expiry date indicated on the box label.

PROTOCOL

Protein preparation and labeling conditions

The labeling process is pH and concentration sensitive. Make sure that the pH of the buffer is 9.0 and that the concentration of the protein is at 16.67 µM (e.g. 2.5 mg/mL for an antibody).

Depending on the molecular weight of the target, each kit enables the labeling of 25 µg to 250 µg (15 kDa to 150 kDa MW respectively) of molecule with an initial molar ratio of 5 d2 per molecule.

Concentration of the molecule to be labeled should be determined at its maximum absorption using the corresponding molar extinction coefficient, i.e. for an antibody:

$$[\text{antibody}] \text{ mole/L} = \frac{\text{OD}_{280\text{nm}}}{210,000}$$

Where 210,000 is the molar extinction coefficient (ϵ) in $\text{cm}^{-1}\text{M}^{-1}$ of a typical IgG at 280nm.

Labeling procedure

The complete procedure is described alongside and takes approximately 1h.

The column supplied with this kit cannot dry out. The volume of eluent recovered at each step corresponds exactly to the volume loaded. Wait until the elution of each step is completed before starting the next one.

Add 100 μL of molecule at 16.67 μM (1.67 nmoles)
(i.e.: 250 μg for an antibody)
directly into the d2 vial



Mix thoroughly with a vortex for 2 min



Incubate for 45 min at RT

The reaction mixture is ready for purification

Equilibrate the column by passing 10 mL of elution buffer
(4 x 2.5 mL) through it
(approx 20 min)

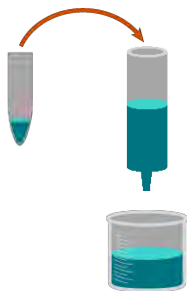
Discard the eluent collected (10 mL)



The column is ready for separation

Conjugate purification:

Add the reaction mixture (100 μL)
onto the column



Discard the eluent
collected (100 μL)

Add 600 μL of elution buffer
onto the column



Discard the eluent
collected (600 μL)

Add 400 μL of elution buffer
onto the column



Collect the 400 μL eluent
(conjugate fraction)

Preparation of the degree of labeling

Measure the absorbance of the purified conjugate in a cuvette with 1cm pathlength at both 280nm (OD_{280nm}) and 650nm (OD_{650nm}).

Determination of the concentration of molecule labeled in the sample (example of an Ab):

$$\text{Molecule concentration (M)} = \frac{OD_{280nm} - (OD_{650nm}/28)}{210,000}$$

- Where 210,000 is the molar extinction coefficient (ϵ) in $\text{cm}^{-1}\text{M}^{-1}$ of a typical IgG at 280nm and 28 a correction factor for the d2's contribution to the absorbance at 280nm.
- The massic (mg/ml) extinction coefficient (ϵ) in $\text{cm}^{-1}\text{M}^{-1}$ of a typical IgG at 280nm is 1.4 and MW is 150,000.

Determination of the final concentration of dye in the sample:

$$\text{Dye concentration} = \frac{OD_{650nm}}{250,000}$$

- Where 250,000 is the molar extinction coefficient (ϵ) in $\text{cm}^{-1}\text{M}^{-1}$ of the d2 dye at 650nm

Determination of the final molecular ratio (MRf) between molecule and dye in the conjugate:

$$\text{Moles of d2 per mole of molecule (MRf)} = \frac{\text{Dye concentration (M)}}{\text{Molecule concentration (M)}}$$

In general, for IgG, the labeling procedure allows 70% of initial material to be recovered and we find that labeling with 2 to 4 moles of d2 per mole of antibody is optimal.

Preparation of the conjugate stock solution

Divide the conjugate into aliquots and freeze at -20°C .

Conjugate Working concentration recommendation

Conjugate working concentration will depend on the assay considered. For a d2-labeled antibody, in 384-well low volume plate format, 20ng of antibody per well (i.e. 0.13 pmol/well) may be used as a starting point. Further assay optimization will require the conjugate titration by assessing a range of antibody concentrations ranging from 1ng to 100 ng per well.

Conjugate storage conditions and handling

Divide the stock solution into suitably sized aliquots and store at -20°C. Avoid repeated thaw/freeze cycles.

Recommendations

- Always store the d2 under a desiccated atmosphere. d2 is stable for 3 months at -60°C or below.
- Strictly follow the instructions. Always start working with a 5-fold amount of d2 per molecule to be labeled to ensure an efficient coupling (labeling efficacy is concentration dependent).
- Impure molecules or material stabilized with carrier proteins cannot be labeled reliably, unless pre-purified.
- Do not elute more than 400 µL of conjugate fraction (step 4).
- Do not re-use the column.

REACH European regulations and compliance

This product and/or some of its components include a Triton concentration of 0.1% or more and as such, it is concerned by the REACH European regulations. We recommend researchers using this product to act in compliance with REACH and in particular: to only use the product for in vitro research in appropriate and controlled premises by qualified researchers, ii) to ensure the collection and the treatment of subsequent waste, and iii) to make sure that the total amount of Triton handled does not exceed 1 ton per year.

This product contains material of biologic origin. Use for research purposes only. Do not use in humans or for diagnostic purposes. The purchaser assumes all risk and responsibility concerning reception, handling and storage.



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF

Manufactured by Cisbio Bioassays - Parc Marcel Boiteux - 30200 Codolet - FRANCE

www.revvity.com

revvity

Revvity, Inc.
940 Winter Street
Waltham, MA 02451 USA

(800) 762-4000
www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2023, Revvity, Inc. All rights reserved.