



Atlas Labeling Kit and Reagents

Product Overview

The Following Kit contains necessary reagents for performing photoinduced proximity labelling using the Atlas system. The procedures are adopted from adapted from Geri et. al., Microenvironment Mapping via Dexter Energy Transfer on Immune Cells, Science (2020). The following kit provides enough antibody conjugate to perform 10 proteomics experiments.

Contents of Atlas Kit

Iridium Photocatalyst (5mM in DMSO) (Red cap – Catalog # IR01), 1 x 8 μ L

Diazirine-PEG3-biotin (Purple cap – Catalog # - DZ01) (MW= 644.7 g/mol), 1 x 10mg

Azido PEG3 NHS ester (Orange cap – Catalog # - CR01), 1 x 18ug

Conjugation Buffer A (DPBS) (Clear cap – Catalog # - BS01), 1 x 5mL

Conjugation Buffer B (1M sodium bicarbonate) (White cap – Catalog # - BS02), 1 x 100 μ L

DMSO (Green cap – Catalog # - OS01), 1 x 100 μ L

Labeling Buffer (DPBS) (Yellow cap - Catalog # - LDSB001), 1 x 15mL

Fluorescent Standard (Catalog # -RF01), 1 x 250 uL



Dexterity Pharma

Additional Materials Required

The following items are not included in the Kit:

Primary antibody

Secondary antibody

Isotype antibody

Milli-Q Water

Vortexer

Centrifuge

LoBind Protein tubes (1.5 mL)

15 mL conical centrifuge tubes

PhotoReactor can be purchased through efficiency aggregators:

<https://www.efficiencyaggregators.com/photoreactor-2/>

<https://www.sigmaaldrich.com/catalog/product/sial/z744035?lang=en®ion=US>

BCA assay (Pierce)

Fluorescence/Absorbance Plate Reader

Important Product Information

- Diazirine-based reagents are both heat and light sensitive. When not in use store in the dark at 4 °C.
- Diazirine-based reagents undergo non-specific labeling when exposed to light <450nm.
- High-intensity blue light exposure should be avoided at all times.
- For best results, Iridium photocatalyst-conjugated antibodies should be prepared and used immediately. Remaining antibody-conjugates can be stored for up to 1 week at 4 °C.
- NHS esters are extremely moisture sensitive and should be equilibrated to room temperature prior to opening.



Dexterity Pharma

- NHS esters are amine reactive and so all amine containing buffers and additives must be removed prior to ligation (i.e. Tris / sodium azide).
- To avoid contamination of samples, always wear gloves, use ultrapure LC/MS grade reagents, and work in a clean work area.

Preparation of Photocatalyst Antibody conjugate

1. Equilibrate azido PEG3 NHS ester (CR01) to room temperature. Add 10 μ L DMSO (OS01) to vial containing azido PEG3 NHS ester (CR01). Vortex to dissolve, centrifuge for 1 minute at 1000xg to collect contents.
2. Add 8 μ L of the DMSO solution of CR01 to IR01 and incubate at room temperature in the dark for 1 h.
3. To a 1.5 mL centrifuge tube, add 300 μ g of your desired antibody, diluted to 150 μ L with Conjugation buffer A.

If antibody is already at 2 mg/mL, there is no need to dilute further with Conjugation buffer A. If antibody is less than 2 mg/mL, the reaction may be less efficient. Antibody solution can be made more concentrated using protein ultracentrifugation concentrators (not included).

4. To the solution of antibody from step 3, add 10 μ L Conjugation Buffer B (1M NaHCO₃, BS02). Mix thoroughly with a pipette. Then, add 17 μ L of the solution from step 2 (i.e CR01 + IR01 solution).
5. Mix thoroughly with pipette and allow the mixture to incubate at 4 °C for 2 h in the dark with end-over-end rotation.
6. Snap off the plastic cap at the bottom of the desalting column (DC01) and loosen the cap at the top of the column. Place the column in an empty 15 mL centrifuge tube. Centrifuge at 1000g for 2 minutes to remove the storage buffer. Discard eluent.
7. Add 1 mL Conjugation Buffer A (BS01) to the top of the column (DC01) from step 5. Centrifuge at 1000g for 2 minutes. Discard eluent.
8. Repeat Step 7.
9. Place desalting column in a new 15 mL centrifuge tube and load the antibody-Ir conjugation reaction onto the column bed. Centrifuge at 1000g for 2 minutes. Collect the eluent.
10. Transfer the eluent to a new Protein Lo-bind tube.



Dexterity Pharma

11. Measure protein concentration using BCA assay (Pierce) / Bradford 595 protein assay (BioRad) according to manufacturer's instructions.
12. Measure fluorescence (385 nm excitation, 495 nm emission) using a plate reader with a black-bottom 384-well plate. Compare against an 8-point calibration curve made from the supplied reference compound (RF01) made using 2-fold serial dilutions in DPBS. The reference compound is 100 μ M of an iridium complex with identical spectroscopic properties as the clickable iridium photocatalyst; fluorescence when compared against the calibration curve will provide the concentration of iridium in the antibody conjugate sample.
13. Divide the concentration of iridium in the antibody conjugate sample by the concentration of the antibody measured through BCA assay (mg/mL divided by the molecular weight of the antibody employed, ca. 150,000 daltons). This number is the average number of iridium photocatalysts per antibody. Typical values range from 2 to 18.
14. Adjust protein concentration to 1 mg/mL using DPBS.



Dexterity Pharma

Photolabeling of cells for quantitative LC/MS/MS analysis

1. Remove cells (recommended 20 million cells per replicate) from culture and pellet according to manufacturer's guidelines (e.g. 300xg for 4 min at 4 °C). From here on, keep cells at 4 °C at all times.
2. Remove supernatant and wash cells with cold DPBS (2 x 1mL) and resuspend in cold DPBS (1mL) containing 25µg targeting primary antibody in a Protein LoBind tube.
3. Incubate cells on a rotisserie for 30 min at 4 °C.
4. Centrifuge cells and remove supernatant.
5. Wash cells twice with 1 mL cold DPBS by gentle resuspension using a pipette followed by centrifugation.
6. Suspend cell pellet gently using a pipette in 1 mL cold DPBS containing 25 µg of photocatalyst conjugated goat α-mouse secondary IgG prepared in step 24.
7. Incubate cells on a rotisserie for 30 min at 4 °C.
8. During this time, add 62 µL of DMSO to vial containing Diazirine-PEG3-biotin (DZ01) and vortex to dissolve. To prepare 250 µM Diazirine-PEG3-biotin solution in DPBS, dilute Diazirine-PEG3-biotin solution to cold DPBS at a dilution of 1:1000 and vortex thoroughly. Keep solution on ice, protected from light.
9. Centrifuge cells and remove supernatant.
10. Wash cells twice with 1 mL cold DPBS by gentle resuspension using a pipette followed by centrifugation.
11. Suspend cell pellet gently using a pipette in 1 mL of cold DPBS containing 250 Diazirine-PEG3-biotin.
12. The samples were placed in the photoreactor (BPR200) and irradiated at 100% intensity (10 minutes in BPR200, 3 minutes in Penn PhD Photoreactor M2).
13. After irradiation, centrifuge cells and remove supernatant.
14. Wash cells twice with 1 mL cold DPBS by gentle resuspension using a pipette followed by centrifugation.

***Cells are now ready for downstream processing, immunoprecipitation, and LC/MS/MS analysis based upon standard procedures and/or user's best practice
For an example, see Oakley et al. Proc. Natl. Acad. Sci. U. S. A. 2022, 119, e2203027119***



Trouble Shooting:

Problem	Possible cause	Solution
Low Iridium concentration on antibody	Primary amine-based buffer used during NHS ligation	Use non-primary amine based buffer (1 M NaHCO ₃ in DPBS)
	Incorrect buffer pH	Ensure pH ~ 8.0–8.5
	Too much sample used	Use 4 µL NHS ester solution
	Click reagents decomposed	Use fresh Click reagents avoiding exposure to air
	Reagents hydrolyzed	Reconstitute NHS ester in anhydrous DMSO
	Antibody decomposition	Reduce amount of NHS ester used for azide ligation
Low labelling efficiency	Diazirine probe decomposition	Store in the dark at –20 °C until use
	Antibody conjugate decomposed	Use antibody conjugate immediately following preparation, avoid storage for extended periods. Storage in 10% glycerol in DPBS if any precipitation occurs.
	Low intensity irradiation	Ensure biophotoreactor is set to 100% intensity during irradiation
	Antibody has low affinity	Use a different primary antibody

Additional Information:

Ask a qualified Proteomics professional for further information regarding data analysis.



Dexterity Pharma

Dexterity Pharma LLC Limited Warranty

Limited Use Label License:

The purchase of this product conveys to the Buyer the non-exclusive, non-transferable right to use the Product and components of the Product for internal research purposes of the Buyer only, regardless of whether the Buyer is an academic or for-profit entity or some other type of entity. This Product is not available for use, and should not be used, in diagnostic procedures. Dexterity Pharma LLC warrants that all products in this Product have passed quality control testing. A sample Product from the same lot was randomly chosen and demonstrated reproducible photoaffinity labelling of a protein target using conjugation to a known antibody to said target. This Product is not intended for human use and Dexterity Pharma LLC is not liable for injuries that may occur from the improper use of this product. Do not ingest or otherwise consume any material included in this Pit. Unless otherwise expressly stated on the Product or in the Documentation accompanying the Product, the Product is intended for research only and is not to be used for any other purpose, including without limitation, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses, or any type of consumption by or application to humans or animals.

By opening and using this Product, the Buyer acknowledges and expressly agrees to be bound by the terms and conditions of this Limited Use Label License and Limited Warranty. This product is warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts (“Documentation”). No claim of suitability for use in applications regulated by FDA is made. This product is not designed to diagnose, treat, cure, or prevent any disease. The one-year Limited Warranty provided herein is valid only when the Product is used by properly trained individuals, and when the Product is subjected to normal, proper and intended usage. Unless otherwise stated in the Documentation, this Limited Warranty is limited to one-year from date of shipment. This warranty does not extend to anyone other than Buyer. The Buyer may not transfer this warranty to any other individual or entity under any circumstances. Any assignment, sale or transfer of this warranty shall immediately terminate all liability of Dexterity Pharma LLC for the Product, and the warranties contained herein. Any model or sample furnished to Buyer is merely illustrative of the general type and quality of goods and does not represent that any Product will conform to such model or sample.

The Buyer must notify Dexterity Pharma LLC of any claims under this Limited Warranty, provide proof of purchase with such notice, and provide the Product thought to be defective back to Dexterity Pharma LLC, within thirty (30) days of the Buyer discovering, or when buyer should have discovered, the Product as nonconforming, at [P.O. Box 353 / 331 N Post Rd /

Princeton Junction, NJ 08550]. Failure of the Buyer to notify Dexterity Pharma LLC, provide proof of purchase, and return such Product to Dexterity Pharma LLC in accordance with this procedure shall constitute a waiver by Buyer of any and all claims of any kind with respect thereto and immediately extinguish all liability of Dexterity Pharma LLC under this warranty and render Buyer’s warranty claim null and void. Any Product returned by Buyer to Dexterity



Dexterity Pharma

Pharma LLC shall be returned to Dexterity Pharma LLC's place of business stated above, freight prepaid, accompanied or preceded by Buyer's particularized statement of the claimed defect in the Product.

TO THE MAXIMUM EXTENT PERMITTED BY APPLICABLE LAW, DEXTERITY PHARMA, LLC, EXPRESSLY DISCLAIMS ANY OTHER WARRANTIES OTHER THAN THE ONE PROVIDED HEREIN, EXPRESS OR IMPLIED, STATUTORY OR OTHERWISE, INCLUDING WITHOUT LIMITATION, IMPLIED WARRANTIES OF MERCHANTABILITY, FITNESS FOR ANY PARTICULAR PURPOSE, TITLE, QUALITY, ACCURACY, QUIET ENJOYMENT AND/OR NON INFRINGEMENT OF THIRD PARTY RIGHTS, AND ANY IMPLIED WARRANTIES ARISING FROM COURSE OF DEALING OR PERFORMANCE. UNDER NO CIRCUMSTANCES WILL DEXTERITY PHARMA LLC BE LIABLE FOR ANY LOSS OR INJURY OR ANY DIRECT, INDIRECT, INCIDENTAL, SPECIAL, RELIANCE, CONSEQUENTIAL, EXEMPLARY OR PUNITIVE DAMAGES WHATSOEVER, INCLUDING, WITHOUT LIMITATION, ANY OTHER LOSS, DAMAGE OR EXPENSE OF ANY KIND FOR INJURY TO PERSON OR PROPERTY, FOR LOSS OF BUSINESS, REVENUE, PROFITS, GOODWILL, BUSINESS INTERRUPTION, LOSS OF BUSINESS INFORMATION, LOSS OF PRIVACY, FAILURE TO MEET ANY DUTY, OR NEGLIGENCE, ARISING OUT OF OR IN ANY WAY RELATED TO THE PRODUCT UNDER ANY LEGAL THEORY INCLUDING WITHOUT LIMITATION BREACH OF CONTRACT, STRICT LIABILITY OR TORT (INCLUDING NEGLIGENCE, OR ANY OTHER ACTION BY ANY PARTY), EVEN IF DEXTERITY PHARMA LLC HAS BEEN ADVISED OF THE POSSIBILITY OF SUCH DAMAGES.

THIS LIMITED WARRANTY IS CONDITIONAL UPON THE REASONABLE OPINION OF DEXTERITY PHARMA LLC, IN ITS SOLE DISCRETION, THAT THERE EXISTS A DEFECT IN THE PRODUCT. IF UPON EXAMINATION BY DEXTERITY PHARMA LLC, THE PRODUCT IS SHOWN TO HAVE A DEFECT UNDER THE ONE-YEAR WARRANTY PERIOD, THE BUYER'S SOLE AND EXCLUSIVE REMEDY IS LIMITED TO REPAIR, REPLACEMENT OF OR REFUND FOR THE NONCONFORMING PRODUCT, TO BE DETERMINED AT SELLER'S SOLE OPTION AND DISCRETION, AND SUCH REMEDY SHALL BE BUYER'S ENTIRE AND EXCLUSIVE REMEDY. DEXTERITY PHARMA LLC'S MAXIMUM LIABILITY SHALL NOT EXCEED THE PURCHASE PRICE OF THE PRODUCT FOUND TO BE DEFECTIVE.

DEXTERITY PHARMA LLC SHALL NOT BE LIABLE UNDER ANY CIRCUMSTANCES FOR, AND THERE IS NO OBLIGATION TO REPAIR, REPLACE OR REFUND FOR THE PRODUCT AS THE RESULT OF:

- (I) ACCIDENT, DISASTER OR EVENT OF FORCE MAJEURE, INCLUDING WITHOUT LIMITATION WAR (WHETHER DECLARED OR NOT); ARMED CONFLICT OR SERIOUS THREAT OF SAME (INCLUDING BUT NOT LIMITED TO HOSTILE ATTACK, BLOCKADE, MILITARY EMBARGO), HOSTILITIES, INVASION, ACT OF A FOREIGN ENEMY, EXTENSIVE MILITARY MOBILIZATION; CIVIL WAR, RIOT, REBELLION, REVOLUTION, MILITARY OR USURPED POWER, INSURRECTION, CIVIL COMMOTION OR DISORDER,



Dexterity Pharma

MOB VIOLENCE, ACT OF CIVIL DISOBEDIENCE; ACT OF GOD OR NATURAL DISASTER (SUCH AS EARTHQUAKE, HURRICANE, LIGHTNING STRIKE, TORNADO, HAIL, ANY OTHER VIOLENT STORM, CYCLONE, TYPHOON, BLIZZARD, VOLCANIC ACTIVITY, LANDSLIDE, TIDAL WAVE, TSUNAMI, FLOOD, DAMAGE OR DESTRUCTION BY LIGHTNING, DROUGHT); ACT OF TERRORISM, SABOTAGE OR PIRACY; PLAGUE, EPIDEMIC, PANDEMIC, OUTBREAKS OF INFECTIOUS DISEASE OR ANY OTHER PUBLIC HEALTH CRISIS, INCLUDING QUARANTINE OR OTHER EMPLOYEE RESTRICTIONS; EXPLOSION, FIRE, DESTRUCTION OF MACHINES, EQUIPMENT, FACTORIES AND OF ANY KIND OF INSTALLATION, PROLONGED BREAK-DOWN OF TRANSPORT, TELECOMMUNICATION OR ELECTRIC CURRENT; GENERAL LABOR DISTURBANCE SUCH AS BUT NOT LIMITED TO BOYCOTT, STRIKE AND LOCK-OUT, GO-SLOW, OCCUPATION OF FACTORIES AND PREMISES; SHORTAGE OR INABILITY TO OBTAIN CRITICAL MATERIAL OR SUPPLIES TO THE EXTENT NOT SUBJECT TO THE REASONABLE CONTROL OF DEXTERITY PHARMA LLC; OR ANY OTHER EVENT OR CONDITION BEYOND THE REASONABLE CONTROL OF DEXTERITY PHARMA LLC;

- (II) MISUSE, IMPROPER USE, ABUSE, FAULT OR NEGLIGENCE OF OR BY BUYER;
- (III) BEING ACCIDENTALLY OR INTENTIONALLY DAMAGED, HAVING NOT BEEN PROPERLY MAINTAINED OR OPERATED, OR HAS BEEN ALTERED OR MODIFIED IN ANY WAY;
- (IV) USE OF THE PRODUCT IN A MANNER FOR WHICH IT WAS NOT DESIGNED;
- (V) IMPROPER SHIPPING, STORAGE OR HANDLING OF THE PRODUCT;
- (VI) DAMAGE TO ANY PERSONAL OR REAL PROPERTY FROM USE OF THE PRODUCT.
- (VII) ANY COSTS, EXPENSES OR DAMAGES, INCLUDING LABOR OR MATERIAL, INCURRED DURING USE OF THE PRODUCT OR PARTS THEREOF; OR
- (VIII) FAILURE OF THE PRODUCT OR PARTS THEREOF FROM EXTERNAL FORCES SUCH AS, BUT NOT LIMITED TO, VANDALISM, VEHICULAR OR OTHER IMPACT.

IN NO EVENT SHALL THE FOREGOING LIMITED WARRANTY EXTEND TO ANY PRODUCT IN ANY WAY CAUSED OR ALLOWED TO BE, OR INSTALLED, OPERATED



Dexterity Pharma

OR USED IN SUCH A MANNER AS TO BE, SUBJECT OR EXPOSED TO CONDITIONS OF MISUSE, ABUSE OR ACCIDENT. ANY ACTION FOR BREACH OF THIS LIMITED WARRANTY OR ANY OTHER ACTION AGAINST DEXTERITY PHARMA LLC RELATING TO OR ARISING OUT OF THE PRODUCT, SHALL BE BROUGHT WITHIN ONE-YEAR AFTER ANY SUCH CAUSE OF ACTION ACCRUES. REPAIR OR REPLACEMENT OF THE PRODUCT UNDER THIS LIMITED WARRANTY SHALL NOT EXTEND THE DURATION OF THE LIMITED WARRANTY WITH RESPECT TO SUCH REPAIRED OR REPLACED PRODUCT BEYOND THE STATED WARRANTY PERIOD.

THIS LIMITED WARRANTY GIVES THE ORIGINAL BUYER OF THE PRODUCT ONLY SPECIFIC LEGAL RIGHTS. YOU MAY ALSO HAVE OTHER RIGHTS WHICH MAY VARY BY EACH JURISDICTION.