

AlphaScreen®

## Conjugation Kit

Product number: 6760000K

Lot Number: 3225063

Product Format: 6760000K 2 x2mg

1mg translates into 2,000 assay points is based on a final bead concentration of 20 µg/mL in a 25 µL/well reaction volume.

Manufacturing date: 11/6/2023

Document version: 1

### Kit Components

Component	6760000K
Unconjugated Acceptor Beads at 20 mg/mL in 0.1M MES, pH 6.0	2 x 50 µL (6762001)
Streptavidin Donor Beads at 5 mg/mL in 25mM Hepes, 0.1M NaCl, 0.05% proclin, pH 7.4	2 x 200 µL (6760002S)

### Product Information

**Stability:** This product is stable for at least 12 months from the manufacturing date if used and stored under recommended conditions.

**Storage:** Store undiluted at 4°C protected from light. Freeze-thaw is not recommended and can cause the beads to form aggregates.

### Conjugation Protocol

**Note:** The following procedure is recommended only as a general guideline for bead conjugation to IgG molecules. It is highly recommended that conditions are further optimized as required for the application for which the beads are intended.

#### Materials:

**Aldehyde beads:** (Provided)

**Sodium cyanoborohydride solution:** Dissolve 25 mg of sodium cyanoborohydride (Sigma #15,615-9 or equivalent) in 1 mL Milli-Q® H<sub>2</sub>O or equivalent. **Note:** Prepare fresh.

**MES buffer (0.1 M pH 6.0):** Dissolve 2.13g MES (2-[N-Morpholino]ethanesulfonic Acid) (Sigma # M-5287 or equivalent) in 80 mL Milli-Q® H<sub>2</sub>O or equivalent. Adjust the pH to 6.0 with 10 N NaOH and make up to 100 mL total volume. Filter through a 0.2 µm filter.

**1% Tween-20:** Dilute 10% Tween-20 (Pierce #28320 or equivalent) 1/10 in MES buffer.

**CMO solution (0.3 M pH 5.0):** Dissolve 65 mg CMO (carboxymethoxylamine hemihydrochloride) (Sigma #C1,340-8 or equivalent) in 1 mL Milli-Q® H<sub>2</sub>O or equivalent. Adjust the pH to 5.0 with 10 N NaOH. **Note:** Prepare fresh.

**TRIS buffer (0.1 M pH 8.0):** Dissolve 1.21g Tris (ICN #819638 or equivalent) in 80mL deionised H<sub>2</sub>O. Adjust the pH to 8.0 with HCl and make up to 100mL total volume. Filter through a 0.2µm filter.

**Conjugation:**

1. To each tube (1 mg beads) provided add:
  - 12.5 µL 1% Tween-20
  - 0.05 mg antibody (This will provide a 20:1 bead:antibody wt:wt ratio. As antibodies are available in different concentrations depending on the supplier, 50 µL of a 1 mg/mL solution is recommended)
  - 10µL Sodium cyanoborohydride solution
2. Make up to a total volume of 200 µL with 0.1 M MES pH 6.0. Incubate in the dark for 48 hours at 37°C.

**Blocking:**

Add 10 µL CMO solution. Incubate in the dark at 37°C for 1 hour.

**Purification:**

1. Add 190 µL 0.1 M Tris buffer pH 8.0.
2. Centrifuge at 13,000 g for 30 min at 4°C.
3. Remove the supernatant with a micropipette.
4. Add 1 mL 0.1 M Tris buffer pH 8.0 to pellet.
5. Vortex and sonicate\*.
6. Centrifuge at 13,000 g for 30 min at 4°C.
7. Repeat steps 3 through 6.
8. Add 200 µL of an appropriate buffer\*\* (5 mg/mL final concentration).
9. Vortex and sonicate\*.
10. Store in the dark at 4°C

*\*Sonicate on ice at 10 pulses of 1 second (0.5 second ON/0.5 second OFF) using a Misonix XL sonicator (Model CL4) and HS419 microtip probe.*

*\*\* Either PBS or 25 mM Hepes, 100 mM NaCl, pH 7.4 is recommended.*

Please visit our website for additional information on AlphaScreen technology at [www.revvy.com](http://www.revvy.com)

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. 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.

6760000K-R Rev 01

revvity