

DBCO–Azide Conjugation

OVERVIEW

Dibenzocyclooctyne (DBCO) reagents enable bioorthogonal “click” reactions with azide-functionalized molecules through Strain-Promoted Alkyne–Azide Cycloaddition (SPAAC). This copper-free reaction forms a stable triazole product under mild conditions, making it ideal for bioconjugation, labeling, and chemical biology applications.

Key Features

- **Biocompatible:** No cytotoxic copper catalyst required (suitable for in vivo applications).
- **Mild Conditions:** Efficient conjugation in aqueous buffer at low or room temperature.
- **Long-Term Stability:** Both DBCO and azide moieties remain stable for extended periods.
- **High Efficiency:** Triazole formation typically proceeds in quantitative yield.
- **Specific & Bioorthogonal:** Azide reacts only with DBCO in the presence of other functional groups (e.g., $-\text{NH}_2$, $-\text{SH}$, $-\text{COOH}$).
- **UV Traceability:** DBCO absorbs strongly at ~ 310 nm, allowing reaction monitoring by UV–Vis.

Notes

- **Solvent Conditions:** Although DBCO–Azide reactions can proceed in partially organic media (e.g., up to 20% DMSO in aqueous buffer), for antibody conjugations we typically recommend standard aqueous buffers (e.g., PBS) without sodium azide, since azide will deplete DBCO.
- **PEG Linkers:** DBCO or azide reagents containing PEG arms offer enhanced hydrophilicity.

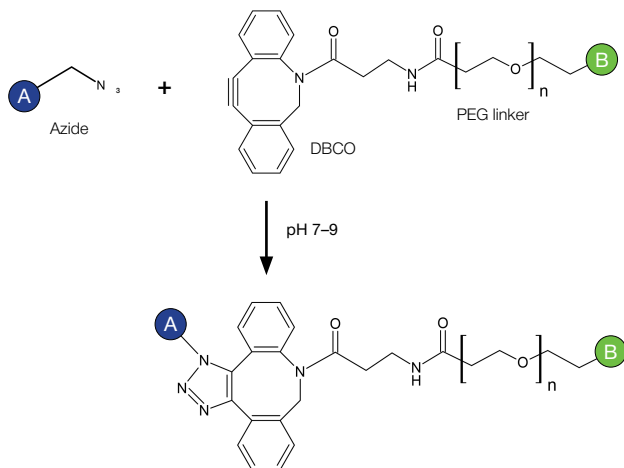


Figure 1. Schematic of SPAAC (Strain-Promoted Azide-Alkyne Cycloaddition) ligation reaction, in which a triazole bond is formed between an azide molecule and DBCO with high efficiency and specificity under mild conditions.

GENERAL LIGATION WORKFLOW

1. Azide Activation of Biomolecule A

- Attach an azide moiety to Biomolecule A.

2. DBCO Activation of Biomolecule B

- Separately, conjugate a DBCO reagent (e.g., DBCO–NHS ester) to Biomolecule B.

3. Click Reaction

- Mix the DBCO-functionalized Biomolecule #1 with the azide-functionalized Biomolecule #2 to form the stable triazole conjugate.

EXAMPLE PROTOCOL: ANTIBODY–OLIGO CONJUGATION

Antibody Activation with DBCO

1. Reaction Setup

- Dissolve DBCO–NHS ester in DMSO at 10 mM.
- Use antibody at 1–10 mg/mL in a suitable buffer (e.g., PBS, pH ~ 7.4). (In many SE-labeling protocols, 2 mg/mL or higher is standard; a range of 1–10 mg/mL is acceptable.)
- Mix the antibody with a 20–30-fold molar excess of DBCO–NHS.
- Keep the final DMSO content at less than 20%.

2. Incubate

- React at room temperature for 60 minutes.

3. (Optional) Quench

- Typically, SE conjugations do not require a specific quench step and proceed directly to purification. However, if desired, add ~ 10 μL of 100 mM glycine (in water) to neutralize unreacted DBCO–NHS.
- Incubate for 15 minutes.

Characterize the DBCO–Antibody (Degree of Substitution)

To calculate the average number of DBCO molecules per antibody (sometimes referred to as DOS or DBCO/Ab), measure the absorbance of the purified conjugate at 280 nm and 309 nm and use the following formula:

$$\text{DBCO} / \text{Ab} = \frac{A_{309} \times \epsilon_{\text{Ab}}}{A_{280} - (\text{CF} \times A_{309}) \times \epsilon_{\text{DBCO}}}$$

Where:

- $\epsilon_{\text{DBCO}} = 12,000 \text{ M}^{-1} \text{ cm}^{-1}$ at 309 nm
- $\epsilon_{\text{Ab}} = 210,000 \text{ M}^{-1} \text{ cm}^{-1}$ at 280 nm (for IgG)
- A_{309} = absorbance at 309 nm
- A_{280} = absorbance at 280 nm
- CF = 0.955 (correction factor for DBCO absorbance bleed at 280 nm)

EXAMPLE PROTOCOL: ANTIBODY-OLIGO CONJUGATION CONT.

Click Chemistry with Azide-Oligonucleotide

1. Set Up Reaction

- Mix the DBCO-antibody with a 2–4-fold molar excess of azide-modified oligonucleotide (or azide-labeled dye).
- Recommended solvent: PBS or similar buffer with up to 20% DMSO. Avoid sodium azide.

2. Incubate

- Incubate 2–4 hours at room temperature (or overnight at 4 °C) to ensure complete reaction.

3. Validation

- Check conjugate formation via SDS-PAGE. The antibody-oligo conjugate should show a higher molecular weight band than the unmodified antibody.

4. Purification

- Remove excess oligonucleotide or dye by liquid chromatography (e.g., reverse-phase or ion-exchange HPLC) or another appropriate method.

EXAMPLE DATA ANALYSIS AND FIGURES

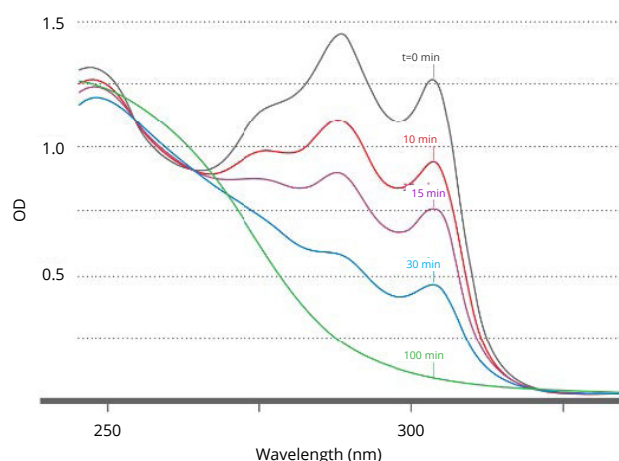


Figure 2. Real-time UV-Vis monitoring of the SPAAC (DBCO-Azide) reaction at various time points (0–100 min). The initial black spectrum (t=0 min) shows a pronounced DBCO peak near 310 nm, which progressively diminishes as the reaction proceeds. By 100 min (green line), the 310 nm peak is significantly reduced, indicating extensive consumption of DBCO.

APPENDIX

Table 1. Ordering information for AAT Bioquest azide and DBCO products

Cat. No.	Product	Quantity
12642	10XHis azide	1 mg
20332	2',3'-cGAMP azide	100 µg
20330	2',3'-cGAMP-DBCO conjugate	100 µg
21243	5-(and-6)-Carboxy SNARF-1 Azide	1 mg
131	5-FAM Azide	10 mg
950	5-FAM Azide	100 mg
486	5-TAMRA azide	5 mg
960	5-TAMRA azide	50 mg
133	6-FAM Azide	10 mg
955	6-FAM Azide	100 mg
240	6-HEX azide	5 mg
248	6-JOE azide	5 mg
217	6-NED azide	1 mg
236	6-OG488 Azide	1 mg
494	6-ROX azide	5 mg
490	6-TAMRA azide	5 mg
965	6-TAMRA azide	50 mg
244	6-TET azide	5 mg
219	6-VIC Azide	1 mg
12628	6XHis azide	1 mg
12623	6XHis-DBCO Conjugate	1 mg
26019	Acridinium Azide	1 mg
508	AMCA Azide	1 mg
70203	ATTO 390 azide	1 mg
70205	ATTO 390 DBCO	1 mg
70213	ATTO 425 azide	1 mg
70215	ATTO 425 DBCO	1 mg
2817	ATTO 488 azide	1 mg
2819	ATTO 488 DBCO	1 mg
2802	ATTO 488 PEG3 azide	1 mg
70223	ATTO 495 azide	1 mg
70225	ATTO 495 DBCO	1 mg
2845	ATTO 514 azide	1 mg
2847	ATTO 514 PEG4 DBCO	1 mg
2824	ATTO 532 PEG3 azide	1 mg
2821	ATTO 532 PEG4 DBCO	1 mg
70233	ATTO 550 azide	1 mg
70235	ATTO 550 DBCO	1 mg
70243	ATTO 590 azide	1 mg
70245	ATTO 590 DBCO	1 mg
2840	ATTO 594 PEG3 azide	1 mg
2842	ATTO 594 PEG4 DBCO	1 mg
70253	ATTO 610 azide	1 mg

APPENDIX CONT.

Cat. No.	Product	Quantity
70255	ATTO 610 DBCO	1 mg
70263	ATTO 620 azide	1 mg
70265	ATTO 620 DBCO	1 mg
70273	ATTO 633 azide	1 mg
70275	ATTO 633 DBCO	1 mg
2832	ATTO 647 PEG4 DBCO	1 mg
2835	ATTO 647N azide	1 mg
2837	ATTO 647N DBCO	1 mg
2806	ATTO 647N PEG3 azide	1 mg
2807	ATTO 647N PEG4 DBCO	1 mg
70283	ATTO 655 PEG3 azide	1 mg
70285	ATTO 655 PEG4 DBCO	1 mg
70293	ATTO 680 PEG3 azide	1 mg
70295	ATTO 680 PEG4 DBCO	1 mg
70303	ATTO 700 PEG3 azide	1 mg
70305	ATTO 700 PEG4 DBCO	1 mg
2412	AzoDye-1 Azide	1 mg
2413	AzoDye-1-PEG10-Azide	1 mg
2432	AzoDye-2 Azide	1 mg
2431	AzoDye-2 C2 Azide	1 mg
2433	AzoDye-2-PEG10 Azide	1 mg
2482	AzoDye-3 Azide	1 mg
2483	AzoDye-3 PEG10 Azide	1 mg
20423	BAPTA Azide	1 mg
20425	BAPTA DBCO	1 mg
12568	BG Azide	1 mg
3023	Biotin Azide	5 mg
3020	Biotin C2 Azide	5 mg
3025	Biotin-PEG2-azide *CAS 945633-30-7*	5 mg
3019	Biotin-PEG3-azide *CAS 875770-34-6*	5 mg
20613	Cal-520® azide	100 µg
6909	CDPI3 Azide [Minor Groove Binder Azide]	1 mg
943	Cy3B azide	1 mg
945	Cy3B DBCO	1 mg
143	Cyanine 3 azide [equivalent to Cy3® azide]	1 mg
970	Cyanine 3 azide [equivalent to Cy3® azide]	25 mg
980	Cyanine 3.5 azide [equivalent to Cy3.5® azide]	1 mg
981	Cyanine 3.5 azide [equivalent to Cy3.5® azide]	25 mg
153	Cyanine 5 azide [equivalent to Cy5® azide]	1 mg
975	Cyanine 5 azide [equivalent to Cy5® azide]	25 mg

Cat. No.	Product	Quantity
178	Cyanine 5.5 azide [equivalent to Cy5.5® azide]	1 mg
163	Cyanine 7 azide [equivalent to Cy7® azide]	1 mg
2010	DABCYL-DBCO	5 mg
4529	DBCO NHS Ester	10 mg
4534	DBCO PEG4 Amine	25 mg
4527	DBCO Sulfo-NHS Ester	5 mg
4528	DBCO Sulfo-NHS Ester	25 mg
920	DBCO-Cy3	1 mg
923	DBCO-Cy5	1 mg
4530	DBCO-PEG4-NHS Ester	1 mg
4526	DBCO-PEG5-NHS Ester	1 mg
3504	Digoxigenin azide	1 mg
72710	FastClick™ 5-FAM Azide	1 mg
72712	FastClick™ 5-TAMRA Azide	1 mg
72711	FastClick™ 6-FAM Azide	1 mg
72714	FastClick™ 6-ROX Azide	1 mg
72713	FastClick™ 6-TAMRA Azide	1 mg
72801	FastClick™ Biotin Azide	1 mg
72700	FastClick™ Cy3 Azide	1 mg
72702	FastClick™ Cy5 Azide	1 mg
72704	FastClick™ Cy7 Azide	1 mg
72800	FastClick™ Digoxigenin (DIG) Azide	1 mg
72730	FastClick™ XFD350 Azide	1 mg
72733	FastClick™ XFD405 Azide	1 mg
72735	FastClick™ XFD488 Azide	1 mg
72737	FastClick™ XFD555 Azide	1 mg
72740	FastClick™ XFD647 Azide	1 mg
72745	FastClick™ XFD750 Azide	1 mg
17616	Helixyte™ Green Azide	1 mg
17674	Hoechst 33342 azide	1 mg
985	ICG azide	1 mg
1094	iFluor® 405 azide	1 mg
1000	iFluor® 488 azide	1 mg
1093	iFluor® 555 azide	1 mg
1095	iFluor® 594 azide	1 mg
1091	iFluor® 647 azide	1 mg
1363	iFluor® 790 Azide	1 mg
20342	IP1 Azide	100 µg
50637	Isotonitazene Azide	1 mg
832	Methylene Blue Azide	1 mg
1690	mFluor™ Violet 450 Azide	1 mg
24215	MycLight™ PMA [Propidium Monoazide]	1 mg
24216	MycLight™ PMA [Propidium Monoazide] *5 mM Water Solution*	100 µL
12605	NTA Azide	1 mg

APPENDIX CONT.

Cat. No.	Product	Quantity
5303	Phalloidin-PEG4-DBCO	100 µg
39057	Psoralen MOP Azide	1 mg
39062	Psoralen TMP Azide	1 mg
3403	ReadiLeave™ Reversible Biotin Azide	1 mg
16883	ReadiUse™ Preactivated Streptavidin Azide	1 mg
21125	Rhod-4™ azide	100 µg
39004	SDA azide	1 mg
50007	Semicarbazide (SEM)-BSA Conjugate	1 mg
50009	Semicarbazide (SEM)-HRP Conjugate	1 mg
50008	Semicarbazide (SEM)-KLH Conjugate	1 mg
484	Texas Red® azide *Single Isomer*	5 mg
2236	Tide Fluor™ 1 azide [TF1 azide]	5 mg
2252	Tide Fluor™ 2 azide [TF2 azide]	1 mg
2254	Tide Fluor™ 3 azide [TF3 azide]	1 mg
2300	Tide Fluor™ 4 azide [TF4 azide]	1 mg
2275	Tide Fluor™ 5WS azide [TF5WS azide]	1 mg
2302	Tide Fluor™ 6WS azide [TF6WS azide]	1 mg
2304	Tide Fluor™ 7WS azide [TF7WS azide]	1 mg
2306	Tide Fluor™ 8WS azide [TF8WS azide] *Near Infrared Emission*	1 mg
2188	Tide Quencher™ 1 azide [TQ1 azide]	5 mg
2211	Tide Quencher™ 2 azide [TQ2 azide]	5 mg
2231	Tide Quencher™ 3 azide [TQ3 azide]	5 mg
2068	Tide Quencher™ 4WS azide [TQ4WS azide]	1 mg
2070	Tide Quencher™ 4WS-DBCO [TQ4WS-DBCO]	1 mg
2088	Tide Quencher™ 5.1WS azide [TQ5.1WS azide]	1 mg
2082	Tide Quencher™ 5WS azide [TQ5WS azide]	1 mg
2097	Tide Quencher™ 6WS azide [TQ6WS azide]	1 mg
2120	Tide Quencher™ 7.1WS azide [TQ7.1WS azide]	1 mg
2128	Tide Quencher™ 7.2WS azide [TQ7.2WS azide]	1 mg
2112	Tide Quencher™ 7WS azide [TQ7WS azide]	1 mg
2136	Tide Quencher™ 8WS azide [TQ8WS azide]	1 mg
1446	trFluor™ Eu DBCO *europium complex*	100 µg
70002	XFD350 azide	1 mg
70004	XFD350 PEG4 DBCO	1 mg
70013	XFD405 azide	1 mg
70015	XFD405 PEG4 DBCO	1 mg
70024	XFD430 azide	1 mg

Cat. No.	Product	Quantity
70026	XFD430 PEG4 DBCO	1 mg
1701	XFD488 azide *Same Structure to Alexa Fluor™ 488 azide*	1 mg
1813	XFD488 PEG4 DBCO	1 mg
70044	XFD514 azide	1 mg
70046	XFD514 PEG4 DBCO	1 mg
1717	XFD532 azide	1 mg
1719	XFD532 PEG4 DBCO	1 mg
70054	XFD546 azide	1 mg
70056	XFD546 PEG4 DBCO	1 mg
1723	XFD555 azide	1 mg
1725	XFD555 PEG4 DBCO	1 mg
70133	XFD568 azide	1 mg
70135	XFD568 PEG4 DBCO	1 mg
1729	XFD594 azide	1 mg
1731	XFD594 PEG4 DBCO	1 mg
70064	XFD610 azide	1 mg
70066	XFD610 PEG4 DBCO	1 mg
70074	XFD633 azide	1 mg
70076	XFD633 PEG4 DBCO	1 mg
70084	XFD635 azide	1 mg
70086	XFD635 PEG4 DBCO	1 mg
1896	XFD647 Azide	1 mg
1734	XFD647 PEG4 DBCO	1 mg
70094	XFD660 azide	1 mg
70096	XFD660 PEG4 DBCO	1 mg
70104	XFD680 azide	1 mg
70106	XFD680 PEG4 DBCO	1 mg
70114	XFD700 azide	1 mg
70116	XFD700 PEG4 DBCO	1 mg
1738	XFD750 azide	1 mg
1740	XFD750 PEG4 DBCO	1 mg
70123	XFD790 azide	1 mg
70125	XFD790 PEG4 DBCO	1 mg

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.