



# Antibody-Oligonucleotide All-in-One™ Conjugation Kit

## A-9202

[Product Images](#)



## Short Description

---

The Antibody-Oligonucleotide All-in-One Conjugation Kit uses patented SoluLINK<sup>®</sup> bioconjugation technology to link an antibody to an oligonucleotide. A succinimidyl-4-formylbenzamide (S-4FB)-modified oligonucleotide and succinimidyl-6-hydrazino-nicotinamide (S-HyNic)-modified antibody form an antibody-oligonucleotide conjugate. This kit is specifically designed to conjugate 100 µg of antibody to any amino-modified oligonucleotide in the 20-80 nucleotide range.

### Key Features:

- Any mammalian monoclonal or polyclonal IgG isotype antibody can be conjugated to the oligonucleotide and purified with approximately 4 hours of total hands-on time.
- SoluLINK bioconjugation technology forms a stable and UV-traceable bis-aryl hydrazone bond with measurable absorbance at 354 nm, which has a molar extinction coefficient of 29,000 L mol<sup>-1</sup> cm<sup>-1</sup>.
- The conjugation reaction is spectrophotometrically quantifiable due to the formation of a chromophoric conjugate bond.

## Additional Information

---

Conjugation Targets	Antibody
Label/Modifier Type	Oligonucleotide
Reactivity	Amine
Recommended Storage	2° - 8°C – Do Not Freeze
Applications	Antibody Labeling, In Situ Proximity Ligation

## Products in this set

---



**Antibody-Oligonucleotide All-in-One™ Conjugation Kit**

