



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

## A-9002

[Product Images](#)



## Short Description

---

The HRP-antibody All-in-One Conjugation Kit contains everything needed to make two HRP-antibody conjugates from 100 µg of any user-supplied antibody. Included in the kit are purification columns to remove unconjugated HRP and antibody, which produces conjugates with superior levels of detection and low non-specific binding.

## Additional Information

---

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

## Products in this set

---



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

