



## Label-Free Biomolecular Interaction Analysis

### Octet® Regeneration Buffer Kit

#### Quickly Screen for Biosensor Regeneration Conditions

- Contains 4 ready-to-use buffers:
  - 10 mM Glycine, pH 1.0, 15 mL
  - 10 mM Glycine, pH 1.7, 15 mL
  - 40 mM Phosphoric Acid, 15 mL
  - 80 mM Phosphoric Acid, 15 mL
  
- Suitable for various Octet® biosensor types and a wide range of ligands and analytes

# Octet® Regeneration Buffer Kit

To determine accurate affinity values, the ligand surface must remain consistent across runs in the same study. This means all analyte molecules must unbind from the ligand prior to assessing subsequent samples or analyte concentrations. When measuring low-affinity interactions (high nM to low  $\mu$ M), it is possible to allow the dissociation phase to continue until all analyte has unbound from the ligand and the buffer baseline response is achieved. This is not a practical approach when measuring high-affinity interactions (pM to low nM), so a chemical intervention is required for regeneration. The Octet® Regeneration Buffer Kit provides a set of ready-to-use buffers for a quick screening of different regeneration conditions for various Octet® biosensor types and a wide range of ligands and analytes. Regeneration allows for biosensor re-use and provides a cost-effective solution for analyzing a large number of samples or ligand-analyte pairs or in sequence.

---

Item	Order Code
------	------------

---

Octet® Regeneration Buffer Kit	18-5171
--------------------------------	---------



Octet® 10mM Glycine pH 1.7, 100 mL (standalone buffer)	18-1184
---	---------



## Germany

Sartorius Lab Instruments GmbH & Co.  
KG Otto-Brenner-Strasse 20  
37079 Goettingen  
Phone +49 551 308 0

## USA

Sartorius Corporation  
565 Johnson Avenue  
Bohemia, NY 11716  
Phone +1 888 OCTET 75  
Or +1 650 322 1360



For further contacts, visit

[www.sartorius.com/octet-support](http://www.sartorius.com/octet-support)

Specifications subject to change without notice.

Copyright Sartorius Lab Instruments GmbH & Co. KG.

For Research Use Only.

Status: 05 | 2024