

## Product datasheet

# Development Enzyme Mix ab154151

### Overview

---

<b>Product name</b>	Development Enzyme Mix
<b>Product overview</b>	Development Enzyme Mix for <a href="#">ab65620</a>

### Properties

---

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	1 units
Development Enzyme Mix (lyophilized)	1 vial

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

### Our Abpromise to you: Quality guaranteed and expert technical support

---

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
  
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

### Terms and conditions

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors