

# NEW » LAB STARTUP PROGRAM

## Save more with each product purchase.

Purchase up to 20,000CHF of Promega products over a 12 month time period and you can save up to 50% in product credits and rewards. Kick off your savings by spending a minimum of 2,500CHF and receive a 700CHF product purchase credit and a 50CHF breakfast basket for your lab.

2,500CHF Spending Level 30% SAVINGS	5,000CHF Spending Level 35% SAVINGS	10,000CHF Spending Level 40% SAVINGS	15,000CHF Spending Level 45% SAVINGS	20,000CHF Spending Level 50% SAVINGS
Purchase 2,500CHF of products for your new lab and receive 750CHF worth of rewards. >>> <b>A breakfast basket valued at 50CHF</b> >>> <b>A credit of 700CHF</b>	Purchase an additional 2,500CHF of products for a total of 5,000CHF and get 1,000CHF of rewards. >>> <b>A breakfast basket valued at 75CHF</b> >>> <b>A credit of 925CHF</b>	Purchase an additional 5,000CHF of products for a total of 10,000CHF and get 2,250CHF of rewards. >>> <b>A breakfast basket valued at 175CHF</b> >>> <b>A credit of 2,075CHF</b>	Additional savings with another purchase 5,000CHF of products, 15,000CHF total, and get 2,750CHF of rewards. >>> <b>A breakfast basket valued at 300CHF</b> >>> <b>A credit of 2,450CHF</b>	Just spend 5,000CHF more on products and get 3,250CHF of rewards. >>> <b>A breakfast basket valued at 400CHF</b> >>> <b>A credit of 2,850CHF</b>
<b>750CHF Total Savings Value</b>	<b>1,750CHF Total Savings Value</b>	<b>4,000CHF Total Savings Value</b>	<b>6,750CHF Total Savings Value</b>	<b>10,000CHF Total Savings Value</b>

Spend 20,000CHF over a 12-month period to get the maximum savings. >>>



## Terms of Program

By submitting this application form, you agree to the following terms and conditions: Lab has opened within the past 12 months. Shipping and dry ice charges are 45.00CHF per shipment (excluding VAT). Promega AG reserves the right to apply changes to the New Lab Start-Up Program at any time without prior notice. The New Lab Setup Program can be combined with other special offers and purchases through our Helix consignment stocks. However, the value in CHF of the discount will be automatically deducted from the coupons offered at the various stages of the program. The duration of the program is 12 months. You authorize a Promega AG representative to contact you in order to process your application. Instruments and Custom Assay Services (CAS) are excluded. Prices and specifications subject to change. **Please allow 2-3 weeks to process your application.**



## NEW » LAB STARTUP PROGRAM

[www.promega.com/start-up](http://www.promega.com/start-up)