

Payment information for the customers of CRL Hungary

Our Dear Customer,

We would like to kindly ask you to use one of the following bank data for your payments:

Beneficiary name: Charles River Laboratories Hungary Kft.

Beneficiary Address: Hungary-8200 Veszprém Szabadságpuszta hrsz 028/1

Currency:	EUR
IBAN:	HU38 1080 0007 1000 0000 1540 7026
Account number:	15407026
Sort Code:	10800007
SWIFT Code:	CITIHUHX
Bank address:	CITIBANK EUROPE PLC 80 Váci Út 1133 Budapest Hungary

Currency:	USD
IBAN:	HU62 1080 0007 2000 0000 1540 7034
Account number:	15407034
Sort Code:	10800007
SWIFT Code:	CITIHUHX
Bank address:	CITIBANK EUROPE PLC 80 Váci út 1133 Budapest Hungary

Our customers shall bear wire-transfer fees charged by both sender and intermediate banks up to the arrival of money to the receiver bank. CRL Hungary shall only bear credit-entry related costs charged by the receiver bank.

Please note that we do not accept payments by check.

In connection with invoices please contact: SOL-accountsreceivable@crl.com

Sincerely,



Ernő Duda
Managing Director