

March 2023



Customer Notification Letter -FAO: Accounts Payable & Procurement Department.

Dear Valued Partner,

CHARLES RIVER LABORATORIES HUNGARY KFT wish to notify you of our new banking relationship with Citibank and provide you updated bank details. Please amend your records to reflect the new bank details per the information below at your earliest convenience, and by no later than 30th April 2023.

Change of Bank details

Our **new** bank details are listed below. We have enclosed verification of these new bank details from Citibank directly with this letter.

Name of Company: CHARLES RIVER LABORATORIES HUNGARY KFT
Company Address: SZABADSAGPUSZTA HRSZ 028/1 VESZPREM 8200 HU
Citibank Branch: Citibank Europe plc Hungarian Branch Office
Swift Code: CITIHUHX

IBAN Number	Currency
HU03108000074000000015408006	HUF
HU27108000075000000015408014	EUR
HU42108000079000000015408049	USD

WHAT DO THESE CHANGES MEAN TO YOU?

Please amend your records to ensure all your future payments are credited to our new account as detailed above. Remittances should be emailed to our AR Team HU-AccountsReceivable@crl.com

We want to assist you to make sure these changes are as seamless as possible for you. If you have any queries in relation to the bank changes, please feel free to contact the same **AR Team** HU-AccountsReceivable@crl.com who will be happy to deal with your query.

Yours sincerely

A handwritten signature in blue ink, appearing to read "János Gerencsér".

János Gerencsér
Director of operations
CHARLES RIVER LABORATORIES HUNGARY KFT

Charles River Laboratories Hungary Kft.

8200 Veszprém, Szabadságpuszta Hrsz. 028/1., Hungary • Tel +36 88 545-300 • Fax +36 88 545-301 • www.criver.com