



08 November 2022

Re: Legal Entity Rationalization

Dear Valued Trading Partner:

The purpose of this letter is to formally notify you that Solvo Biotechnológiai Zrt. ("SOLVO") is entering into a Plan of Merger (the "Merger Agreement") with Charles River Laboratories Hungary ("CRL HUNGARY") which shall become effective as of **01 December 2022**. Upon completion of the Merger, SOLVO will cease to exist as a legal entity, and CRL HUNGARY will be the surviving entity.

Being a valued trading partner, your company will notice these changes in all documentation received from SOLVO as of the effective date above. To facilitate these changes in your financial systems, please:

Change the details of Solvo Biotechnológiai Zrt. to the following from the effective date:

Entity Name:	Charles River Laboratories Hungary Kft.
Street:	Szabadságpuszta hrsz 028/1
City:	Veszprém
ZIP code:	8200
Country:	Hungary
EU VAT Number:	HU10268856

Please update your master data file accordingly to reflect these changes.

This change reflects the change of legal entity. The existing facilities of SOLVO in Budapest and Szeged, the website www.solvobiotech.com as well as the SOLVO brand name will remain in use for our products and services.

Should you have any questions related to this notification, please reach out to your primary contact at SOLVO, the SOLVO accounts receivable team by email: SOL-accountsreceivable@crl.com or the SOLVO general inquiry email address: solvobiotech@crl.com.

Sincerely,

Ernő Duda
Managing Director

SOLVO Biotechnológiai Zrt.
6724 Szeged, Szilánk köz 3.
Adószám: 12706978-2-06
(2)