

To our Horizon Discovery Customers, our emails will be changing to Revvity, but our billing information will continue to be Horizon Discovery Biosciences Limited. All email addresses will change to the **@revvity.com** domain, and the first part of the email addresses may also vary per user. We have included some FAQs to help address some of the most common questions, but of course, please reach out if you have any additional questions.

Why did you change your email to @revvity.com and is this for everyone at Horizon Discovery?

As Horizon Discovery is now part of Revvity Inc., we're transitioning to the Revvity email system. This change applies to all Horizon employees as well as all our customer care so all email addresses will switch to the @revvity.com domain.

Do I just change @horizondiscovery.com to @revvity.com in my contact information for Horizon Discovery?

Please note that while you will need to update the domain to @revvity.com, the first part of the email addresses might also change per user, based on the new Revvity email structure. Your Horizon Discovery representative's updated email address will be included in their email signature. During the transition period, we will continue to receive emails sent to both our Horizon Discovery and Revvity addresses.

How do I contact the Customer Care/ Customer Service, Technical Support, or Sales teams? Will the email addresses remain the same?

During the transition period, we will continue to receive emails sent to both our Horizon Discovery and Revvity addresses for a period of time, the following outlines the new emails we ask you to make note of.

	Previously Used Email	New Revvity Email
Horizon D365 Lasernet	D365Laseret@horizondiscovery.com	D365Laseret.horizon@revvity.com
Horizon Discovery Orders	orders@horizondiscovery.com	orders.horizon@revvity.com
Horizon Discovery Sales	sales@horizondiscovery.com	sales.horizon@revvity.com
Horizon Discovery Science Support	technical@horizondiscovery.com	technical.horizon@revvity.com
Horizon Discovery EU Science Support	ts.dharmacon.eu@horizondiscovery.com	ts.dharmancon.eu@revvity.com

Please use our [Contact Us](#) page to send a message directly to our teams.

Do I need to change anything in my ordering or payment information for Horizon Discovery going forward?

No, you don't need to change anything at this time. Your ordering and payment information will remain the same.

Will I still order from the [Horizondiscovery.com](https://horizondiscovery.com/) website?

Yes, for now, please continue to visit <https://horizondiscovery.com/> to place orders, view updates, and access product information. We will notify you when this web address changes.

Do I still use the same phone numbers?

Yes, you can continue to use the same phone numbers. You can expect the same excellent Service from our Customer Care team.

Is anything else changing that I should be aware of?

No, not at this time. We'll make sure to communicate any future company changes to you, but of course if you have any questions, please reach out to our customer care team.

Are product offerings changing?

No, all product offerings will remain the same.