



Temporary Suspension of Reverse Phase Protein Array (RPPA) Services

Dear Customers,

Thank you for using our RPPA Services.

The glass slides we use to create protein arrays from customer supplied lysates have unfortunately been discontinued by our supplier, forcing us to temporarily suspend our service using this format. Over the past few months, we researched alternative suppliers and attempted to optimize experimental conditions to allow use of a new source and avoid interruption of the service. To our regret, this was not possible, as we have not found an acceptable replacement. As soon as we are able to resume our RPPA Services, we will post the announcement to this website.

We deeply value our relationship with our customers and sincerely apologize for any difficulty that may be caused by this development.

In the meantime, if you have any questions at all, please do not hesitate to contact us at info@carnabio.com.

Sincerely,

Carna Biosciences, Inc.