



Product Notice

Date: March 18, 2016

Catalog No. 31400-1

Description: Anti-Human Interferon Beta, Sheep Serum (PAb)

This notice is to inform you that this product has been discontinued due to exhaustion of the supply of this reagent. PBL recommends product 31401-1 Anti-Human Interferon Beta, Sheep Serum (PAb) as a suitable replacement.

<https://www.pblsaysci.com/content/anti-human-interferon-beta-sheep-serum>

Product 31401-1 is offered at the same price as the discontinued product and is suitable for the same applications (tested and suggested).

The principal difference between these two polyclonal antibodies is in the antigen from which they were derived. The replacement polyclonal antibody is derived from sheep immunized with **mammalian cell-expressed** human IFN beta whereas the original product was derived from sheep immunized with *E.coli*-expressed human IFN beta protein.

We thank you for your patience as PBL continues to improve the quality of its products. Please feel free to contact us at techservice@pblsaysci.com if you have any product related questions.