

**TGF-(BETA)1 LAP-D**

**(R58) (LAP**

**Degradates C-**

**Terminus side cut end**

**R58) monoclonal**

Ordering Information

[Order Online »](#)

**antibody**

CAC-CAC-RIK-MA-R58	100µg
--------------------	-------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      -20°C

**Regulatory Status** RUO - Research Use Only

## Product Details

**Application**                      IHC

**Clone**                                18F9-16

**Host**                                 Mouse

**Isotype**                            IgG3Kappa

**Species Reactivity**            Mouse



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)