



A Geno Technology, Inc. (USA) brand name

Immobilized Ficin

For the Generation of Fab & Fc Fragments from Mouse IgG₁

(Cat. # 786-793)



INTRODUCTION	3
ITEM(S) SUPPLIED	
STORAGE CONDITIONS	
IMPORTANT INFORMATION	
ADDITIONAL COMPONENT(S) REQUIRED	
PREPARATION BEFORE USE	
PROCEDURE	5
RELATED PRODUCTS	6

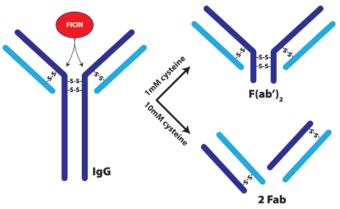
INTRODUCTION

Ficin (or Ficain) (~25,000Da) is a cysteine protease enzyme (EC 3.4.22.3) isolated from fig latex is that has the endopeptidase activity to cleave immunoglobulin G molecules in the hinge reason. Ficin has an effective range of pH4-9.5 with an optimal pH of 6.5 and cleaves bonds that involve uncharged or aromatic amino acids.

Ficin is typically used to cleave mouse IgG_1 as this are difficult to cleave with papain and pepsin. In the presence of 1mM or 10mM cysteine, ficin generates $F(ab')_2$ and $F(ab')_2$ fragments respectively. The Fab and $F(ab')_2$ fragments can be separated from whole IgG and $F(ab')_2$ with either Immobilized Protein A (Cat. # 786-283) or ion exchange chromatography.

Immobilized Ficin is a convenient reagent for producing Fab and $F(ab')_2$ fragments as it avoids the need to remove the ficin enzyme after digestion.

Supplied as a 30% slurry in 50% glycerol, 10mM sodium tetrathionate, 2mM EDTA, pH7.0.



ITEM(S) SUPPLIED

(-) -		
Cat. #	Description	Size
786-793	Immobilized Ficin	5ml resin

STORAGE CONDITIONS

Shipped at ambient temperature. Upon receipt store at 4°C, do NOT freeze.

IMPORTANT INFORMATION

• Activity: 1-1.5mg ficin/ml of resin

Support: 6% Cross-linked Agarose

ADDITIONAL COMPONENT(S) REQUIRED

- Cysteine.HCl
- EDTA
- Sample Buffer: 0.1M Citrate buffer, pH6.0
- Purified, lyophilized IgG or ≥20mg/ml IgG solution
- Wash Buffer: 10mM Tris.HCl, pH7.5

PREPARATION BEFORE USE

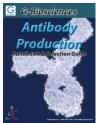
- 10X F(ab')₂ Digestion Buffer: Immediately prior to digestion, add 18.6mg EDTA to 1ml 0.1M Citrate buffer, pH6.0 and once dissolved add 1.8mg Cysteine.HCl to prepare the 10X concentrated buffer.
- 10X Fab Digestion Buffer: Immediately prior to digestion, add 18.6mg EDTA to 1ml 0.1M Citrate buffer, pH6.0 and once dissolved add 18mg Cysteine.HCl to prepare the 10X concentrated buffer.
- 3. *Ficin Activation Buffer:* Dilute the appropriate 10X Digestion Buffer in 0.1M Citrate buffer, pH6.0. Add 1ml Digestion Buffer to 9ml 0.1M Citrate buffer, pH6.0.
- Antibody Preparation: If using an IgG solution, dialyze against the Sample Buffer and concentrate to ~0.5-10mg/ml. Add 100μl appropriate Digestion Buffer to each ml of antibody
 - **NOTE:** We recommend using Tube-O-DIALYZER $^{\text{TM}}$ (Cat. # 786-610 to 786-624) for dialysis to ensure no loss of antibody.
- 5a *Resin Preparation (Column Digestion):* Suspend the resin by gently shaking and inverting the resin. Transfer 2-4ml of the slurry (1-2ml resin) to a suitable column with a wide bore pipette tip. Equilibrate the resin with the addition of 20ml Ficin Activation Buffer. Allow the Ficin Activation Buffer to pass through the column by gravity flow.
- 5b Resin Preparation (Suspension Digestion): Suspend the resin by gently shaking and inverting the resin. Transfer 2-4ml of the slurry (1-2ml resin) to a 50ml tube o with a wide bore pipette tip. Equilibrate the resin with the addition of 20ml Ficin Activation Buffer. Centrifuge at 1,000g for 2-5minutes to pellet the resin, remove the Digestion Buffer.

PROCEDURE

- Add the 1.1ml IgG sample to the activated Immobilized Ficin. Add an additional 0.25ml Activation Buffer to the column to ensure sample fully enters the resin. Seal the tube/column and incubate at 37°C in a high speed shaking waterbath for the indicated time:
 - a. For Mouse IgG₁ Fab fragments incubate for 3-5 hours
 - b. For Mouse IgG_1 F(ab')₂ fragments incubate for 20 hours when using 0.5-3mg/ml antibody
 - For Mouse IgG₁ F(ab')₂ fragments incubate for 40 hours when using 3-10mg/ml antibody
- Centrifuge and collect the fragment containing supernatant or the column flowthrough.
- 3. To separate the Fab fragments from the Fc fragments, use Immobilized Protein A (Cat. # 786-283) or ion exchange. Do not use Protein G as Fab fragments, as well as Fc fragments have some affinity for Protein G.

RELATED PRODUCTS

Download our Antibody Purification Handbook.



 $\underline{\text{http://info.gbiosciences.com/complete-Antibody-Production-handbook}}$

For other related products, visit our website at $\underline{www.GBiosciences.com}$ or contact us.

Last saved: 2/27/2019 CMH

This page is intentionally left blank



www.GBiosciences.com