



130PR

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

FOCUS™ Extraction Buffers

Trial Kit

(Cat. # 786-234)



think proteins! think G-Biosciences www.GBiosciences.com

INTRODUCTION

The FOCUS Extraction Buffers are proprietary modifications of well-studied protein solubilization and IPG-strip re-hydration buffers designed to produce better spot resolution for 2D gel analysis. Depending on the nature of your samples, one of the following buffers will be suitable for most of your applications. The Extraction Buffer-I and II are suitable for most applications; however, for stronger solubilization effects, the Extraction Buffer-III, -IV, -V or -VI may be used. The FOCUS™ Extraction Buffer Trial Kit is designed for those who prefer to experiment for establishing a suitable extraction and solubilization buffer for their protein samples. The trial kit consists of one each of the FOCUS Extraction Buffers I-VI with their corresponding DILUENT.

ITEM(S) SUPPLIED (Cat. # 786-234)

Description	Composition	Size
FOCUS™ Extraction Buffer-I	Urea plus NP-40	For 10ml (5.0g)
FOCUS™ Extraction Buffer-II	Urea plus CHAPS	For 10ml (5.0g)
FOCUS™ Extraction Buffer-III	Urea, Thiourea, ASB-16 plus CHAPS	For 10ml (5.0g)
FOCUS™ Extraction Buffer-IV	Urea, Thiourea, SB 3-10, plus CHAPS	For 10ml (5.0g)
FOCUS™ Extraction Buffer-V	Urea, Thiourea, plus CHAPS	For 10ml (5.0g)
FOCUS™ Extraction Buffer-VI	Urea, Thiourea, NDSB 201, plus CHAPS	For 10ml (5.0g)
DILUENT-I	NP-40	7ml
DILUENT-II	CHAPS	14ml
DILUENT-III	CHAPS	12ml

STORAGE CONDITION

The kit is shipped at ambient temperature. Upon arrival, store FOCUS Extraction Buffers at room temperature and DILUENTS at 4°C.

PROTOCOL

1. Hydrate the supplied FOCUS Extraction Buffers I to VI by adding 5.57ml or 5.0ml of specific DILUENT. Add the specified DILUENT (as per the table on page 2) directly into the corresponding FOCUS Extraction Buffer bottle.
OPTIONAL: Smaller volume can also be prepared by using 1g of dry powder mix with 1ml or 1.15ml specific DILUENT as per table on page 2.
2. Add needed agents such as reducing agents, inhibitors, carrier ampholyte, Bromophenol blue dye, etc.
3. Mix periodically by vortexing and incubate at room temperature until you have a clear solution.

NOTE: Extraction Buffer-IV might require warming for a short period at 25-30°C (water bath).

- Use of only freshly hydrated FOCUS Extraction Buffer is highly recommended. However, any unused buffer can be stored in small aliquots (1-2ml) at -20°C.
- The following table shows the type of DILUENT to use with FOCUS Extraction Buffers I to VI:

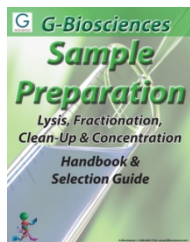
FOCUS™ Extraction Buffer	Volume of DILUENT	DILUENT	W/V Ratio of FOCUS™ Extraction Buffer & DILUENT
FOCUS™ Extraction Buffer-I	5.75ml	I	1g/1.15ml DILUENT-I
FOCUS™ Extraction Buffer-II	5.75ml	II	1g/1.15ml DILUENT-II
FOCUS™ Extraction Buffer-III	5.0ml	III	1g/1ml DILUENT-III
FOCUS™ Extraction Buffer-IV	5.0ml	III	1g/1ml DILUENT-III
FOCUS™ Extraction Buffer-V	5.75ml	II	1g/1.15ml DILUENT-II
FOCUS™ Extraction Buffer-VI	5.0ml	III	1g/1ml DILUENT-III

ORDERING INFORMATION

Cat #	Description	Size
786-220	FOCUS™ Extraction Buffer-I (25g) with DILUENT-I (30ml)	For 50ml
786-221	FOCUS™ Extraction Buffer-II (25g) with DILUENT-II (30ml)	For 50ml
786-222	FOCUS™ Extraction Buffer-III (25g) with DILUENT-III (30ml)	For 50ml
786-223	FOCUS™ Extraction Buffer-IV (25g) with DILUENT-III (30ml)	For 50ml
786-219	FOCUS™ Extraction Buffer-V (25g) with DILUENT-II (30ml)	For 50ml
786-233	FOCUS™ Extraction Buffer-VI (25g) with DILUENT-III (30ml)	For 50ml

RELATED PRODUCTS

Download our Sample Preparation Handbook.



<http://info.gbiosciences.com/complete-protein-sample-preparation-handbook/>

For other related products, visit our website at www.GBiosciences.com or contact us.



www.GBiosciences.com