

# BRFF-HPC1™

BRFF-HPC1™ is a complete serum-free medium designed for establishing new cell lines from human prostate tissue [1]. Epithelial cell line, from both benign prostatic hyperplasia and prostatic carcinoma have been established and maintained in this medium. This medium contains dihydrotestosterone. For optimum cell attachment, use FNC Coating Mix®.

## Ordering Information

[Order Online »](#)

AES-0403	500ml
----------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Use/Stability** 4-6 weeks at +4°C

**Short Term Storage** +4°C

**Long Term Storage** -80°C

**Shipping** Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Formulation** Liquid; Ready-to-Use.

**Technical Info / Product Notes** **pH:** 7.2 -7.5 (5% CO<sub>2</sub> atmosphere is recommended for culture)  
**Osmolality:** 270 -300mOsM



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)