



CURRENT CONTAINER LABEL



Toxic. R25 R35



8°C = 30°C

IVD

SYNCHRON® Systems

TBIL

300

100 mL


REF 442745 (Kit)

LOT

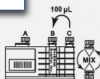
A999999

YYYY-MM-DD

Area Reserved for Instrument Barcode




Beckman Coulter Ireland Inc. 353 91 774068
©2012 Beckman Coulter, Inc.
Beckman Coulter, Inc.,
Brea, CA 92821 U.S.A.



100 µL


B21996 AA




**BECKMAN
Coulter**
Made in U.S.A.

81198123

UPDATED CONTAINER LABEL



DANGER
H401 H302 H314 H317



8°C = 30°C

IVD

SYNCHRON Systems

TBIL

300

100 mL

REF 442745 (Kit)


Country of Origin

YYYY-MM-DD

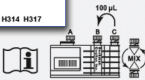
LOT

A999999

Area Reserved for Instrument Barcode




Beckman Coulter Ireland Inc. 353 91 774068
©2014 Beckman Coulter, Inc.
Beckman Coulter, Inc.,
Brea, CA 92821 U.S.A.



100 µL

B21996 XX



**BECKMAN
Coulter**

81198123


CURRENT KIT LABEL

TBIL 2 × 300

[REF] 442745


[LOT] A999999

[EXP] YYYY-MM-DD



+49208697061Q*

+49531170308495000Q*



Toxic. R25 R35 S45 S26 Sodium Nitrite < 10%.

Total Bilirubin Reagent

For use on SYNCHRON Systems for the quantitative determination of total bilirubin in serum or plasma. Contents: 2 × 100 mL Total Bilirubin Reagent. Reactive Ingredients: Sodium Nitrite 1.05 mmol/L, Sodium Acetate 553 mmol/L, Sodium Benzoate 372 mmol/L, Caffeine 181 mmol/L, Sulphanilic Acid 27.8 mmol/L, Hydrochloric Acid 85.4 mmol/L. DO NOT FREEZE.

Reactivo Bilirrubina Total para Sistemas SYNCHRON®. Para uso en los sistemas SYNCHRON para la determinación cuantitativa de bilirrubina total en suero o plasma. Contenido: 2 × 100 mL Reactivo Bilirrubina Total. NO CONGELAR.

Total Bilirubin Reaktifii için SYNCHRON® Sistemleri. SYNCHRON Sistemlerinde serum veya plazmada total bilirubinin kantitatif tayininde kullanım içindir. İçindekiler: 2 × 100 mL Total Bilirubin Reaktifi. DONDURMAYIN.

总胆红素检测试剂 适用于 SYNCHRON®系统。组成：2 × 100 mL 总胆红素检测试剂。不可冷冻。生产商：贝克曼库尔特有限公司，美国加利福尼亚州，Brea市，S. Kraemer大道250号，邮编92821 电话：(001)714-955-5321


[IVD] **CE**

[REF] 442745


[EC REP] Beckman Coulter Ireland, Inc.
Mervyn Business Park,
Mervyn, Galway, Ireland 303 91 774050

© 2014 Beckman Coulter, Inc.
Beckman Coulter, Inc.
250 S. Kraemer Blvd., Brea, CA 92821 U.S.A.

≈ 8°C ≈ 30°C



[REF] 442745

 techdocs.beckmancoulter.com

[EC REP] Beckman Coulter Ireland, Inc.
Mervyn Business Park,
Mervyn, Galway, Ireland 303 91 774050

© 2014 Beckman Coulter, Inc.
Beckman Coulter, Inc.
250 S. Kraemer Blvd., Brea, CA 92821 U.S.A.

**BECKMAN
COULTER**
MADE IN U.S.A.

V 091328


UPDATED KIT LABEL

TBIL 2 × 300

[REF] 442745

[LOT] YYYY-MM-DD

[EXP] A999999



+49208697061Q*

+49531170308495000Q*

[IVD]

DANGER

14691 H302 H314 H317 Sulphanilic Acid 0.1 - 1%, Sodium Acetate 1 - 10%, Sodium Benzoate 1 - 10%, Sodium Nitrite 1 - 10%, Hydrochloric Acid 0.1 - 1%.

Total Bilirubin Reagent

For use on SYNCHRON Systems for the quantitative determination of total bilirubin in serum or plasma. Contents: 2 × 100 mL Total Bilirubin Reagent. Reactive Ingredients: Sodium Nitrite 1.05 mmol/L, Sodium Acetate 653 mmol/L, Sodium Benzoate 372 mmol/L, Caffeine 181 mmol/L, Sulphanilic Acid 27.8 mmol/L, Hydrochloric Acid 85.4 mmol/L. DO NOT FREEZE.

Reactivo Bilirrubina Total para Sistemas SYNCHRON®. Para uso en los sistemas SYNCHRON para la determinación cuantitativa de bilirrubina total en suero o plasma. Contenido: 2 × 100 mL Reactivo Bilirrubina Total. NO CONGELAR.

Total Bilirubin Reaktifii için SYNCHRON® Sistemleri. SYNCHRON Sistemlerinde serum veya plazmada total bilirubinin kantitatif tayininde kullanım içindir. İçindekiler: 2 × 100 mL Total Bilirubin Reaktifi. DONDURMAYIN.

生产商：贝克曼库尔特有限公司，美国加利福尼亚州，Brea市，S. Kraemer大道250号，邮编92821 电话：(001)714-955-5321

[IVD] **CE**


[REF] 442745

[SDS]


[EC REP] Beckman Coulter Ireland, Inc.
Mervyn Business Park,
Mervyn, Galway, Ireland 303 91 774050

© 2014 Beckman Coulter, Inc.
Beckman Coulter, Inc.
250 S. Kraemer Blvd., Brea, CA 92821 USA

≈ 8°C ≈ 30°C



[REF] 442745

 techdocs.beckmancoulter.com

[EC REP] Beckman Coulter Ireland, Inc.
Mervyn Business Park,
Mervyn, Galway, Ireland 303 91 774050

© 2014 Beckman Coulter, Inc.
Beckman Coulter, Inc.
250 S. Kraemer Blvd., Brea, CA 92821 USA

**BECKMAN
COULTER**

06 091328

(81) 15898590217969
(17) 168401
(10) 00328



MADE IN USA