

GEBRUIK VAN DRI[®] ETHYLGLUCURONIDE BECKMAN COULTER DxC 500 AU[®]



Beckman Coulter-reagens REF E0750310

Deze toepassing is bedoeld voor de kwalitatieve en semikwantitatieve bepaling van ethylglucuronide in menselijke urine tot een grenswaarde van 500 ng/ml.



Uitsluitend voor diagnostisch gebruik in vitro
Alleen op voorschrift

Beoogd gebruik



De informatie in dit toepassingsblad is bedoeld als aanvulling op de bijsluiter. Raadpleeg de bijsluiter voor informatie over beoogd gebruik, opslag van reagentia, reagensvoorbereiding, monsterverzameling, monstervoorbereiding, opslag van monsters, kwaliteitscontrole en aanvullende prestatiegegevens.

Bestelinformatie

| Artikel | Grootte | Bestelnummer Beckman Coulter |
|--|------------|------------------------------|
| DRI-ethylglucuronideanalyse | 1 x 18 ml | E0750310 |
| Negatieve kalibrator voor DRI-ethylglucuronide | 1 x 25 ml | E0750311 |
| Kalibrator tot 100 voor DRI-ethylglucuronide | 1 x 10 ml | E0750312 |
| Kalibrator tot 500 voor DRI-ethylglucuronide | 1 x 10 ml | E0750313 |
| Kalibrator tot 1000 voor DRI-ethylglucuronide | 1 x 10 ml | E0750314 |
| Kalibrator tot 2000 voor DRI-ethylglucuronide | 1 x 10 ml | E0750315 |
| Controlemiddel 375 voor DRI-ethylglucuronide | 1 x 25 ml | E0750316 |
| Controlemiddel 625 voor DRI-ethylglucuronide | 1 x 25 ml | E0750317 |
| AU-fles | 20 x 30 ml | 63094 |

Technische ondersteuning

Neem voor technische ondersteuning contact op met uw plaatselijke Beckman Coulter-vertegenwoordiger.

Vervolg op de volgende pagina

Reagensopslag Raadpleeg de bijsluiters voor informatie over de opslag van reagentia.

Analysatorprocedure Raadpleeg de gebruikershandleidingen voor informatie over de werking van de analyser. Raadpleeg de bijsluiters voor de volledige bereiding van het reagens.

Laat het reagens 15 minuten in gekoelde toestand (2 tot 8 °C) stabiliseren voordat u het in AU-flessen giet. Voeg R1-reagens en R2-reagens toe aan de juiste AU-flessen, zoals weergegeven in de onderstaande tabel:

| | AU-reagensfles | |
|-------------------------------------|------------------|------------------|
| DRI-set met ethylglucuronideanalyse | R1-compartiment | R2-compartiment |
| Reagens R1, antistof/substraat | Eén fles (30 ml) | |
| Reagens R2, enzymconjugaat | | Eén fles (30 ml) |

Waarschuwing: deze reagentia moeten op vaste posities worden geprogrammeerd. Gebruik de Thermo-reagensflessen niet rechtstreeks op de AU-analyser.

Interpretatie van resultaten en gegevens Resultaten voor monsters worden afgedrukt in ng/ml.

Vorbereiding van het monster Raadpleeg de bijsluiters voor de volledige preparatie van het monster. De bijsluiters vindt u op de website van Thermo Fisher Scientific:
www.thermofisher.com

Kalibratie Gebruik de DRI-ethylglucuronidekalibrators. De kalibrators zijn vloeibaar en klaar voor gebruik. Raadpleeg de bijsluiters voor de concentratie van elke kalibrator.

Naam reagens: DRI-ethylglucuronideanalyse (alleen kwalitatief tot 500 ng/ml)
REF E0750310 DxC 500 AU urine-instellingen
Naam kalibrator: DRI-ethylglucuronidekalibrator REF E0750313

Reagens-ID 558

TEST CONFIGURATION & CHEMISTRY DETAILS

| | | | | |
|---------------------------------|--------------------------------------|---|--|--|
| Assay Name | Test | Rev | Discipline | Chemistry |
| Test ID | <input type="text" value="ETG500"/> | | Calculated Result | <input type="checkbox"/> |
| LIS Code | <input type="text" value="ETG500"/> | | Result Type | <input type="text" value="Qualitative"/> |
| UNITS AND RANGE SETTINGS | | | | |
| Use Settings from | <input type="text" value="None"/> | Units | <input type="text" value="None"/> | Decimal Places |
| | | | | <input type="text" value="x"/> |
| Test Kind | <input type="text" value="General"/> | Revision | <input type="text" value="01"/> | <input checked="" type="checkbox"/> Multi Reagent Switch |
| Reagent Name | <input type="text" value="ETG"/> | Reagent ID | <input type="text" value="558"/> | <input type="checkbox"/> FSE Test |
| ABB Name | <input type="text" value="ETG1N"/> | Parameter Long Name | <input type="text" value="Ethyl Glucur 500 (Q) E0750310 ETG1N Urine"/> | |
| Region | <input type="checkbox"/> US | <input checked="" type="checkbox"/> OUS | <input checked="" type="checkbox"/> AP | <input type="checkbox"/> JP |
| | | | <input checked="" type="checkbox"/> EU | <input type="checkbox"/> Other |

GENERAL PARAMETERS

| | | | | | |
|------------------------|---|---|-----------------------------------|---|---|
| SAMPLE VOLUME | Sample Volume <input type="text" value="25.0"/> μL | Dilution <input type="text" value="0"/> μL | REACTION OD LIMIT | Low <input type="text" value="-2.0000"/> | High <input type="text" value="3.0000"/> |
| | Predilution Rate <input type="text" value="1"/> | | REACTION BLANK OD LIMIT | First: Low <input type="text" value="-2.0000"/> | High <input type="text" value="3.0000"/> |
| REAGENT VOLUME | R1-1 <input type="text" value="57"/> μL | Dilution <input type="text" value="0"/> μL | | Last: Low <input type="text" value="-2.0000"/> | High <input type="text" value="3.0000"/> |
| | R2-1 <input type="text" value="57"/> μL | Dilution <input type="text" value="0"/> μL | ANALYTICAL MEASURING RANGE | Low <input type="text" value="100.00"/> | High <input type="text" value="2000.00"/> |
| WAVELENGTH | Primary <input type="text" value="340"/> nm | Secondary <input type="text" value="410"/> nm | MANUFACTURER FACTOR | A <input type="text" value="1"/> | B <input type="text" value="0"/> |
| METHOD | <input type="text" value="FIXED"/> | | REAGENT ONBOARD STABILITY | <input type="text" value="31"/> Days | <input type="text" value="0"/> Hours |
| REACTION SLOPE | <input type="text" value="+"/> | | LIH INFLUENCE CHECK | <input type="checkbox"/> Perform LIH check | |
| MEASURING POINT | Point 1: First <input type="text" value="13"/> | Last <input type="text" value="17"/> | Lipemia | <input type="text" value="+"/> | |
| | Point 2: First <input type="text"/> | Last <input type="text"/> | Icterus | <input type="text" value="+"/> | |
| Linearity Limit | <input type="text"/> % | | Hemolysis | <input type="text" value="+"/> | |
| Lag Time Check | <input type="checkbox"/> Perform Lag Time Check | | | | |

Naam reagens: DRI-ethylglucuronideanalyse (alleen kwalitatief tot 500 ng/ml)
 REF E0750310 DxC 500 AU urine-instellingen
 Naam kalibrator: DRI-ethylglucuronidekalibrator REF E0750313, vervolg

Reagens-ID 558

CALIBRATION PARAMETERS

| Base Unit | Decimal Place | Unit 1 | Factor 1 | Unit 2 | Factor 2 | Unit 3 | Factor 3 | Unit 4 | Factor 4 |
|-----------|---------------|--------|----------|--------|----------|--------|----------|--------|----------|
| None | 0 | None | 0 | None | 0 | None | 0 | None | 0 |

CALIBRATOR SPECIFIC

Calibration Type

Counts

CALIBRATION OD AND CONCENTRATION PARAMETERS

Use highest calibrator for Upper AMR

Formula

MB Factor

| | Calibrator Name | Conc | Factor Range Low | Factor Range High |
|---------|-----------------|------|------------------|-------------------|
| Point 1 | ETG CAL-3 | 500 | -99999 | 99999 |
| Point 2 | | | | |
| Point 3 | | | | |
| Point 4 | | | | |
| Point 5 | | | | |
| Point 6 | | | | |
| Point 7 | | | | |

Calibrator Name

Positive Cutoff

SLOPE CHECK

Number of Levels

Slope Check

STABILITY AND INTERVAL

Reagent Blank Stability Days Hours

Interval

Calibration Stability Days Hours

Interval

OD DELTA CHECK

Reagent Blank

Calibration

PROZONE CHECK PARAMETERS

Logic Check 1

Check Points
 Point 1
 Point 2
 Point 3

Decision Values
 Value 1
 Value 2
 Value 3

Logic Check 2

Check Points
 Point 1
 Interval

Limit Points
 Limit 1
 Limit 2

Logic Check 3

Check Points
 Point 1
 Interval

Limit Points
 Limit 1
 Limit 2

Decision Values
 Value 1
 Value 2

Check Pattern
 Pattern

Naam reagens: DRI-ethylglucuronidessay (alleen semi-kwantitatief tot 500 ng/ml) REF E0750310 Dx C 500 AU urine-instellingen

Reagens-ID 558

Naam kalibrator: DRI-ethylglucuronidekalibrators REF E0750311, E0750312, E0750313, E0750314, E0750315

TEST CONFIGURATION & CHEMISTRY DETAILS

| | | | | | | |
|--------------------------|-----------------------------|---|--|--|--|--------------------------------|
| Assay Name | Test | Rev | Discipline | Chemistry | | |
| Test ID | ETG500- | | Calculated Result | <input type="checkbox"/> | | |
| LIS Code | ETG500- | | Result Type | Semiquantitative ▼ | | |
| UNITS AND RANGE SETTINGS | | | | | | |
| Use Settings from | None ▼ | Units | ng/mL ▼ | Decimal Places | x.xx ▼ | Urine |
| Test Kind | General ▼ | Revision | 01 | <input checked="" type="checkbox"/> Multi Reagent Switch | | |
| Reagent Name | ETG | Reagent ID | 558 | <input type="checkbox"/> FSE Test | | |
| ABB Name | ETG2N | Parameter Long Name | Ethyl Glucur S/Q E0750310 ETG2N Urine | | | |
| Region | <input type="checkbox"/> US | <input checked="" type="checkbox"/> OUS | <input checked="" type="checkbox"/> AP | <input type="checkbox"/> JP | <input checked="" type="checkbox"/> EU | <input type="checkbox"/> Other |

GENERAL PARAMETERS

| | | | | | | | | | | |
|-----------------|--|------------------|-----------|---|--------|----------------------------|--|---------|------|---------|
| SAMPLE VOLUME | | Sample Volume | 25.0 μL | Dilution | 0 ▼ μL | REACTION OD LIMIT | Low | -2.0000 | High | 3.0000 |
| REAGENT VOLUME | | Predilution Rate | 1 ▼ | | | REACTION BLANK OD LIMIT | First: Low | -2.0000 | High | 3.0000 |
| | | R1-1 | 57 μL | Dilution | 0 μL | | Last: Low | -2.0000 | High | 3.0000 |
| | | R2-1 | 57 μL | Dilution | 0 μL | ANALYTICAL MEASURING RANGE | Low | 100.00 | High | 2000.00 |
| WAVELENGTH | | Primary | 340 nm | Secondary | 410 nm | MANUFACTURER FACTOR | A | 1 | B | 0 |
| METHOD | | | FIXED 1 ▼ | | | REAGENT ONBOARD STABILITY | | 31 Days | | 0 Hours |
| REACTION SLOPE | | | + | | | LIH INFLUENCE CHECK | <input type="checkbox"/> Perform LIH check | | | |
| MEASURING POINT | | Point 1: First | 13 | Last | 17 | Lipemia | + | ▼ | | |
| | | Point 2: First | | Last | | Icterus | + | ▼ | | |
| Linearity Limit | | | | | | Hemolysis | + | ▼ | | |
| Lag Time Check | | | | <input type="checkbox"/> Perform Lag Time Check | | | | | | |

Naam reagens: DRI-ethylglucuronideanalyse (alleen semi-kwantitatief tot 500 ng/ml) REF E0750310 DxC 500 AU urine-instellingen

Reagens-ID 558

Naam kalibrator: DRI-ethylglucuronidekalibrators REF E0750311, E0750312, E0750313, E0750314, E0750315, *vervolg*

CALIBRATION PARAMETERS

| Base Unit | Decimal Place | Unit 1 | Factor 1 | Unit 2 | Factor 2 | Unit 3 | Factor 3 | Unit 4 | Factor 4 |
|-----------|---------------|--------|----------|--------|----------|--------|----------|--------|----------|
| ng/mL | 2 | None | 0 | None | 0 | None | 0 | None | 0 |

CALIBRATOR SPECIFIC

Calibration Type

Counts

Formula

MB Factor

Calibrator Name

Positive Cutoff

SLOPE CHECK

Number of Levels

Slope Check

STABILITY AND INTERVAL

Reagent Blank Stability Days Hours

Interval

Calibration Stability Days Hours

Interval

CALIBRATION OD AND CONCENTRATION PARAMETERS

Use highest calibrator for Upper AMR

| | Calibrator Name | Conc | OD Range Low | OD Range High |
|---------|-----------------|---------|--------------|---------------|
| Point 1 | ETG CAL-1 | 0.00 | -2.00 | 3.00 |
| Point 2 | ETG CAL-2 | 100.00 | -2.00 | 3.00 |
| Point 3 | ETG CAL-3 | 500.00 | -2.00 | 3.00 |
| Point 4 | ETG CAL-4 | 1000.00 | -2.00 | 3.00 |
| Point 5 | ETG CAL-5 | 2000.00 | -2.00 | 3.00 |
| Point 6 | | | | |
| Point 7 | | | | |

OD DELTA CHECK

Reagent Blank
 Calibration

PROZONE CHECK PARAMETERS

Logic Check 1

Check Points
 Point 1
 Point 2
 Point 3

Decision Values
 Value 1
 Value 2
 Value 3

Logic Check 2

Check Points
 Point 1
 Interval

Limit Points
 Limit 1
 Limit 2

Logic Check 3

Check Points
 Point 1
 Interval

Limit Points
 Limit 1
 Limit 2

Decision Values
 Value 1
 Value 2

Check Pattern
 Pattern

Aanvullende informatie

Belangrijk

Omdat Beckman Coulter het reagens niet produceert of kwaliteitscontroles of andere tests uitvoert op afzonderlijke partijen, is Beckman Coulter niet verantwoordelijk voor de kwaliteit van de verkregen gegevens die worden veroorzaakt door de prestaties van het reagens, variaties in reagenspartijen of protocolwijzigingen door de fabrikant.

Transportschade

Neem contact op met uw Beckman Coulter Technical Support Center als dit product bij ontvangst beschadigd is.

© 2023 Thermo Fisher Scientific Inc. Alle rechten voorbehouden. Alle handelsmerken zijn het eigendom van Thermo Fisher Scientific en haar dochterondernemingen, tenzij anders vermeld. AU Series Systems zijn handelsmerken van Beckman Coulter.



B.R.A.H.M.S GmbH, Neuendorfstrasse 25, 16761, Hennigsdorf, Duitsland

Einde