

ANWENDUNG FÜR DRI® OXYCODONE BECKMAN COULTER DxC 500 AU / 500i



Beckman Coulter-Reagenz REF 100248

Diese Anwendung ist für die qualitative und semiquantitative Bestimmung von Oxycodon in humanem Urin vorgesehen.



Nur zur In-vitro-Diagnostik
Verschreibungspflichtig

Verwendungs- zweck



Die Informationen in diesem Anwendungsblatt sind eine Ergänzung zur Packungsbeilage. Die Packungsbeilage enthält Informationen zu Verwendungszweck, Reagenzienlagerung, Aufbereitung des Reagenz, Probensammlung, Probenvorbereitung, Probenlagerung, Qualitätskontrolle sowie weitere Leistungsdaten.

Bestellinformationen

Artikel	Größe	Beckman Coulter- Nachbestellnummer
DRI Oxycodone Assay	1 x 68 ml	100248
DRI Negativkalibrator	1 x 10 ml	1664
DRI Oxycodone 100 ng/ml Kalibrator	1 x 10 ml	100250
DRI Oxycodone 300 ng/ml Kalibrator	1 x 10 ml	100251
DRI Oxycodone 500 ng/ml Kalibrator	1 x 10 ml	100252
DRI Oxycodone 1000 ng/ml Kalibrator	1 x 10 ml	100253
DRI Oxycodone 100 ng/ml Kontrolle	1 x 10 ml	100254
DRI Oxycodone 300 ng/ml Kontrolle	1 x 10 ml	100255
AU Bottle	20 x 30 ml	63094

Technische Unterstützung

Wenden Sie sich für technische Unterstützung an die Beckman Coulter-Vertretung in Ihrer Region.

Fortsetzung auf der nächsten Seite

Reagenzien- lagerung

Informationen zur Reagenzienlagerung sind in der Packungsbeilage zu finden.

Analyse- verfahren

Informationen zum Betrieb des Analysegeräts sind in dessen Gebrauchsanweisung zu finden. Ausführliche Informationen zur Reagenzienvorbereitung sind in der Packungsbeilage zu finden.

Das Reagenz vor dem Füllen in die AU-Fläschchen 15 Minuten bei Kühltemperatur (2 bis 8 °C) äquilibrieren lassen. Reagenz R1 und Reagenz R2 wie in der folgenden Tabelle gezeigt in die entsprechenden AU-Fläschchen geben:

	AU-Reagenzflasche	
DRI Oxycodone Assay-Kit	R1-Fach	R2-Fach
Antikörper-/Substratreagenz R1	1 Fläschchen (30 ml)	
Enzymkonjugatreagenz R2		1 Fläschchen (30 ml)

Achtung: Diese Reagenzien müssen auf festen Positionen programmiert werden. Die Thermo Reagenzflaschen nicht direkt im AU-Analysegerät verwenden.

Ergebnisse und Dateninterpretation

Die Ergebnisse werden in ng/ml ausgegeben.

Probenvorbereitung

Ausführliche Informationen zur Probenvorbereitung sind in der Packungsbeilage zu finden. Die Packungsbeilage ist auf der Website von Thermo Fisher Scientific zu finden:

www.thermofisher.com

Kalibrierung

Verwenden Sie DRI Oxycodon-Kalibratoren. Die Kalibratoren sind flüssig und gebrauchsfertig. Die Konzentration der Kalibratoren ist der Packungsbeilage zu entnehmen.

Name des Reagens DRI Oxycodone Assay (Qualitativ – 100 ng/ml nur Cutoff)
REF 100248 DxC 500 AU / 500i Urine Settings Name des Kalibrators DRI Oxycodone
Kalibrator REF 100250

Reagenz-ID 557

TEST CONFIGURATION & CHEMISTRY DETAILS

Assay Name	Test	Rev	Discipline	Chemistry
Test ID	OXY100		Calculated Result	☐
LIS Code	OXY100		Result Type	Qualitative ▼

UNITS AND RANGE SETTINGS

Use Settings from	None ▼	Units	None ▼	Decimal Places	x ▼	Urine
Test Kind	General ▼	Revision	01	☒ Multi Reagent Switch		
Reagent Name	OXY	Reagent ID	557	☐ FSE Test		
ABB Name	OXY1G	Parameter Long Name	Oxycodone 100 (Q) 100248 OXY1G Urine			

Region ☒US ☒OUS ☒AP ☐JP ☒EU ☐Other

GENERAL PARAMETERS

SAMPLE VOLUME	Sample Volume	14.0 µL	Dilution	0 ▼ µL	REACTION OD LIMIT	Low	-2.0000	High	3.0000
	Predilution Rate	1 ▼			REACTION BLANK OD LIMIT	First: Low	-2.0000	High	3.0000
REAGENT VOLUME	R1-1	87 µL	Dilution	0 µL		Last: Low	-2.0000	High	3.0000
	R2-1	87 µL	Dilution	0 µL	ANALYTICAL MEASURING RANGE	Low	0.00	High	1000.00
WAVELENGTH	Primary	340 nm	Secondary	410 nm	MANUFACTURER FACTOR	A	1	B	0
METHOD	FIXED ▼				REAGENT ONBOARD STABILITY		31 Days		0 Hours
REACTION SLOPE	+				LIH INFLUENCE CHECK	☐ Perform LIH check			
MEASURING POINT	Point 1: First	17	Last	21	Lipemia	+ ▼			
	Point 2: First		Last		Icterus	+ ▼			
					Hemolysis	+ ▼			
Linearity Limit			%						
Lag Time Check			☐ Perform Lag Time Check						

Name des Reagens DRI Oxycodone Assay (Qualitativ – 100 ng/ml nur Cutoff)
REF 100248 Dx C 500 AU / 500i Urine Settings Name des Kalibrators DRI Oxycodone
Kalibrator REF 100250, Fortsetzung

Reagenz-ID 557

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 ☐ Use highest calibrator for Upper AMR

Formula MB Factor

Calibrator Name Positive Cutoff

☐ SLOPE CHECK Number of Levels

Slope Check

CALIBRATION OD AND CONCENTRATION PARAMETERS

	Calibrator Name	Conc	Factor Range Low	Factor Range High
Point 1	OXY CAL-2	100	-999999	99999
Point 2				
Point 3				
Point 4				
Point 5				
Point 6				
Point 7				

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	0
Point 2	0
Point 3	0

 Limit Points

Limit 1	0
Limit 2	27

 Check Pattern

Pattern	Pattern 1
---------	--

☐ Logic Check 2

Decision Values

Value 1	0
Value 2	0
Value 3	0

 Limit Points

Limit 1	0
Limit 2	27

☐ Logic Check 3

Decision Values

Value 1	0
Value 2	0

 Limit Points

Limit 1	0
Limit 2	27

Name des Reagens DRI Oxycodone Assay (Qualitativ – 300 ng/ml nur Cutoff)
REF 100248 DxC 500 AU / 500i Urine Settings Name des Kalibrators DRI Oxycodone
Kalibrator REF 100251

Reagenz-ID 557

TEST CONFIGURATION & CHEMISTRY DETAILS

Assay Name	Test	Rev	Discipline	Chemistry
Test ID	OXY300		Calculated Result	☐
LIS Code	OXY300		Result Type	Qualitative ▼

UNITS AND RANGE SETTINGS

Use Settings from	None ▼	Units	None ▼	Decimal Places	x ▼	Urine
Test Kind	General ▼	Revision	01	☒ Multi Reagent Switch		
Reagent Name	OXY	Reagent ID	557	☐ FSE Test		
ABB Name	OXY2G	Parameter Long Name	Oxycodone 300 (Q) 100248 OXY2G Urine			

Region ☒US ☒OUS ☒AP ☐JP ☒EU ☐Other

GENERAL PARAMETERS

SAMPLE VOLUME	Sample Volume	10.5 μL	Dilution	0 ▼ μL	REACTION OD LIMIT	Low	-2.0000	High	3.0000
	Predilution Rate	1 ▼			REACTION BLANK OD LIMIT	First: Low	-2.0000	High	3.0000
REAGENT VOLUME	R1-1	87 μL	Dilution	0 μL		Last: Low	-2.0000	High	3.0000
	R2-1	87 μL	Dilution	0 μL	ANALYTICAL MEASURING RANGE	Low	0.00	High	1000.00
WAVELENGTH	Primary	340 nm	Secondary	410 nm	MANUFACTURER FACTOR	A	1	B	0
METHOD	FIXED ▼				REAGENT ONBOARD STABILITY		31 Days		0 Hours
REACTION SLOPE	+				LIH INFLUENCE CHECK	☐ Perform LIH check			
MEASURING POINT	Point 1: First	17	Last	21	Lipemia	+	▼		
	Point 2: First		Last		Icterus	+	▼		
					Hemolysis	+	▼		
Linearity Limit		%							
Lag Time Check	☐ Perform Lag Time Check								

Name des Reagens DRI Oxycodone Assay (Qualitativ – 300 ng/ml nur Cutoff)
REF 100248 Dx C 500 AU / 500i Urine Settings Name des Kalibrators DRI Oxycodone
Kalibrator REF 100251, Fortsetzung

Reagenz-ID 557

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

Formula

MB Factor

Calibrator Name

Positive Cutoff

Add

☐ 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	Factor Range Low	Factor Range High
Point 1	OXY CAL-3	300	-999999	99999
Point 2				
Point 3				
Point 4				
Point 5				
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

Limit Points

Limit 1
 Limit 2

Check Pattern

Pattern

☐ 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

Name des Reagens DRI Oxycodone Assay (Semiquantitativ – 100 ng/ml nur Cutoff)

Reagenz-ID 557

REF 100248 DxC 500 AU / 500i Urine Settings

Name des Kalibrators DRI Oxycodone Kalibratoren REF 1664, 100250, 100251, 100252, 100253

TEST CONFIGURATION & CHEMISTRY DETAILS

Assay Name	Test	Rev	Discipline	Chemistry		
Test ID	OXY100-		Calculated Result	☐		
LIS Code	OXY100-		Result Type	Semiquantitative ▼		
UNITS AND RANGE SETTINGS						
Use Settings from	None ▼	Units	ng/mL ▼	Decimal Places	x.xx ▼	Urine
Test Kind	General ▼	Revision	01	☒ Multi Reagent Switch		
Reagent Name	OXY	Reagent ID	557	☐ FSE Test		
ABB Name	OXY3G	Parameter Long Name	Oxycodone 100 S/Q 100248 OXY3G Urine			
Region	☒ US	☒ OUS	☒ AP	☐ JP	☒ EU	☐ Other

GENERAL PARAMETERS

SAMPLE VOLUME	Sample Volume	10.5	µL	Dilution	0	µL	REACTION OD LIMIT	Low	-2.0000	High	3.0000	
	Predilution Rate	1	▼				REACTION BLANK OD LIMIT	First: Low	-2.0000	High	3.0000	
REAGENT VOLUME	R1-1	87	µL	Dilution	0	µL		Last: Low	-2.0000	High	3.0000	
	R2-1	87	µL	Dilution	0	µL	ANALYTICAL MEASURING RANGE	Low	0.00	High	1000.00	
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	☐ Perform LIH check				
MEASURING POINT	Point 1: First	17		Last	21		Lipemia	+	▼			
	Point 2: First			Last			Icterus	+	▼			
							Hemolysis	+	▼			
Linearity Limit			%									
Lag Time Check	☐ Perform Lag Time Check											

Name des Reagens DRI Oxycodone Assay (Semiquantitativ – 100 ng/ml nur Cutoff)

Reagenz-ID 557

REF 100248 DxC 500 AU / 500i Urine Settings

Name des Kalibrators DRI Oxycodone Kalibratoren REF 1664, 100250, 100251, 100252, 100253 Fortsetzung

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 5AB

Counts 2

Formula POLYGONAL

MB Factor

Calibrator Name

Positive Cutoff 100

☒ SLOPE CHECK

Number of Levels 5

Slope Check +

STABILITY AND INTERVAL

Reagent Blank Stability Days

Hours

Interval Bottle

Calibration Stability Days

Hours

Interval Bottle

CALIBRATION OD AND CONCENTRATION PARAMETERS

☐ Use highest calibrator for Upper AMR

	Calibrator Name	Conc	OD Range Low	OD Range High
Point 1	OXY CAL-1	0.00	-2.00	3.00
Point 2	OXY CAL-2	100.00	-2.00	3.00
Point 3	OXY CAL-3	300.00	-2.00	3.00
Point 4	OXY CAL-4	500.00	-2.00	3.00
Point 5	OXY CAL-5	1000.00	-2.00	3.00
Point 6				
Point 7				

OD DELTA CHECK

☐ Reagent Blank☐ Calibration

PROZONE CHECK PARAMETERS

☐ Logic Check 1

Check Points

Point 1	0
Point 2	0
Point 3	0

Decision Values

Value 1	0
Value 2	0
Value 3	0

☐ Logic Check 2

Check Points

Point 1	0
Interval	1

Decision Values

Value 1	0
Value 2	0

☐ Logic Check 3

Check Points

Point 1	0
Interval	1

Decision Values

Value 1	0
Value 2	0

Limit Points

Limit 1	0
Limit 2	27

Limit Points

Limit 1	0
Limit 2	27

Limit Points

Limit 1	0
Limit 2	27

Check Pattern

Pattern Pattern 1

Name des Reagens DRI Oxycodone Assay (Semiquantitativ – 300 ng/ml nur Cutoff)

Reagenz-ID 557

REF 100248 DxC 500 AU / 500i Urine Settings

Name des Kalibrators: DRI Oxycodone Kalibratoren REF 1664, 100250, 100251, 100252, 100253

TEST CONFIGURATION & CHEMISTRY DETAILS

Assay Name	Test	Rev	Discipline	Chemistry		
Test ID	OXY300-		Calculated Result	☐		
LIS Code	OXY300-		Result Type	Semiquantitative ▼		
UNITS AND RANGE SETTINGS						
Use Settings from	None ▼	Units	ng/mL ▼	Decimal Places	x.xx ▼	Urine
Test Kind	General ▼	Revision	01	☒ Multi Reagent Switch		
Reagent Name	OXY	Reagent ID	557	☐ FSE Test		
ABB Name	OXY4G	Parameter Long Name	Oxycodone 300 S/Q 100248 OXY4G Urine			
Region	☒ US	☒ OUS	☒ AP	☐ JP	☒ EU	☐ Other

GENERAL PARAMETERS

SAMPLE VOLUME		REACTION OD LIMIT					
Sample Volume	10.5 µL	Low	-2.0000	High	3.0000		
Predilution Rate	1 ▼						
REAGENT VOLUME		REACTION BLANK OD LIMIT					
R1-1	87 µL	First: Low	-2.0000	High	3.0000		
R2-1	87 µL	Last: Low	-2.0000	High	3.0000		
Dilution		ANALYTICAL MEASURING RANGE					
0 µL		Low	0.00	High	1000.00		
WAVELENGTH		MANUFACTURER FACTOR					
Primary	340 nm	Secondary	410 nm	A	1	B	0
METHOD		REAGENT ONBOARD STABILITY					
FIXED 1 ▼		31 Days				0 Hours	
REACTION SLOPE		LIH INFLUENCE CHECK					
+		☐ Perform LIH check					
MEASURING POINT		Lipemia				+	▼
Point 1: First	17	Last	21	Icterus	+	▼	
Point 2: First		Last		Hemolysis	+	▼	
Linearity Limit						%	
Lag Time Check		☐ Perform Lag Time Check					

Name des Reagens DRI Oxycodone Assay (Semiquantitativ – 300 ng/ml nur Cutoff)

Reagenz-ID 557

REF 100248 Dx C 500 AU / 500i Urine Settings

Name des Kalibrators: DRI Oxycodone Kalibratoren REF 1664, 100250, 100251, 100252, 100253 Fortsetzung

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 5AB

Counts 2

CALIBRATION OD AND CONCENTRATION PARAMETERS

☐ Use highest calibrator for Upper AMR

Formula POLYGONAL

MB Factor

Calibrator Name

Positive Cutoff 300

Add OXY

☒ SLOPE CHECK

Number of Levels 5

Slope Check +

STABILITY AND INTERVAL

Reagent Blank Stability

Days

Hours

Interval Bottle

Calibration Stability

Days

Hours

Interval Bottle

OD DELTA CHECK

☐ Reagent Blank☐ Calibration

	Calibrator Name	Conc	OD Range Low	OD Range High
Point 1	OXY CAL-1	0.00	-2.00	3.00
Point 2	OXY CAL-2	100.00	-2.00	3.00
Point 3	OXY CAL-3	300.00	-2.00	3.00
Point 4	OXY CAL-4	500.00	-2.00	3.00
Point 5	OXY CAL-5	1000.00	-2.00	3.00
Point 6				
Point 7				

PROZONE CHECK PARAMETERS

☐ Logic Check 1

Check Points

Point 1	0
Point 2	0
Point 3	0

Decision Values

Value 1	0
Value 2	0
Value 3	0

☐ Logic Check 2

Check Points

Point 1	0
Interval	1

Decision Values

Value 1	0
Value 2	0

☐ Logic Check 3

Check Points

Point 1	0
Interval	1

Decision Values

Value 1	0
Value 2	0

Limit Points

Limit 1	0
Limit 2	27

Limit Points

Limit 1	0
Limit 2	27

Limit Points

Limit 1	0
Limit 2	27

Check Pattern

Pattern Pattern 1

Weitere Informationen

Wichtiger Hinweis

Da Beckman Coulter weder das Reagenz herstellt noch Qualitätskontrollen oder andere Prüfungen mit einzelnen Chargen durchführt, schließt Beckman Coulter jegliche Haftung für Qualitätsminderungen der gewonnenen Daten aus, die durch die Leistung des Reagenz, Abweichungen zwischen den Reagenzienchargen oder Protokoll-änderungen des Herstellers entstehen.

Versandschaden

Bitte benachrichtigen Sie umgehend das Technical Support Center von Beckman Coulter, wenn Sie dieses Produkt beschädigt erhalten haben.

© 2024 Thermo Fisher Scientific Inc. Alle Rechte vorbehalten. Alle Marken sind Eigentum von Thermo Fisher Scientific und ihren Tochtergesellschaften, falls nicht anders angegeben. Die Systeme der AU-Serie sind eingetragene Marken von Beckman Coulter.



  B.R.A.H.M.S GmbH, Neuendorfstraße 25, 16761, Hennigsdorf, Deutschland

Ende