

APPLICAZIONE DI DRI[®] OXYCODONE BECKMAN COULTER DxC 500 AU / 500i



Reagente Beckman Coulter REF 100248

La presente applicazione è destinata alla determinazione qualitativa e semiquantitativa dell'ossicodone nelle urine umane.



Solo per uso diagnostico in vitro
Solo su prescrizione medica

Uso previsto



Le informazioni fornite nel presente foglio applicativo costituiscono un'integrazione al foglio illustrativo della confezione. Consultare il foglio illustrativo per informazioni sull'uso previsto, la conservazione e preparazione del reagente, il prelievo, la preparazione e la conservazione dei campioni, il controllo di qualità e ulteriori dati sulle prestazioni.

Informazioni per l'ordine

Articolo	Formato	Numero di riordine Beckman Coulter
DRI Oxycodone Assay	1 x 68 mL	100248
DRI Negative Calibrator	1 x 10 mL	1664
DRI Oxycodone 100 ng/mL Calibrator	1 x 10 mL	100250
DRI Oxycodone 300 ng/mL Calibrator	1 x 10 mL	100251
DRI Oxycodone 500 ng/mL Calibrator	1 x 10 mL	100252
DRI Oxycodone 1000 ng/mL Calibrator	1 x 10 mL	100253
DRI Oxycodone 100 ng/mL Control	1 x 10 mL	100254
DRI Oxycodone 300 ng/mL Control	1 x 10 mL	100255
Flacone in AU	20 x 30 mL	63094

Supporto tecnico

Per il supporto tecnico, contattare il rappresentante locale Beckman Coulter.

Continua alla pagina seguente

**Conservazione
dei reagenti**

Per informazioni sulla conservazione dei reagenti, fare riferimento al foglietto illustrativo.

**Funzionamento
dell'analizzatore**

Per informazioni sul funzionamento dell'analizzatore, fare riferimento ai manuali dell'operatore. Per informazioni sulla preparazione completa dei reagenti, fare riferimento al foglietto illustrativo.

Prima di versare in flaconi AU, lasciare equilibrare il reagente per 15 minuti a temperatura refrigerata (2-8 °C). Dispensare il reagente R1 e il reagente R2 in appositi flaconi AU come indicato nella tabella seguente:

Kit DRI Oxycodone Assay	Flacone per reagente AU	
	Scomparto R1	Scomparto R2
Reagente anticorpo/substrato R1	Un flacone (30 mL)	
Reagente enzima-coniugato R2		Un flacone (30 mL)

Avvertenza: questi reagenti devono essere programmati in posizioni fisse. Non utilizzare i flaconi di reagente Thermo direttamente sull'analizzatore AU.

**Risultati e
interpretazione
dei dati**

I risultati verranno stampati in ng/mL.

**Preparazione
dei campioni**

Per informazioni sulla preparazione completa del campione, fare riferimento al foglietto illustrativo. L'inserito del prodotto è disponibile sul sito Web di Thermo Fisher Scientific:

www.thermofisher.com

Calibrazione

Utilizzare i calibratori per ossicodone DRI. I calibratori sono liquidi e pronti all'uso. Per la concentrazione di ciascun calibratore, fare riferimento al foglietto illustrativo.

Nome reagente: DRI Oxycodone Assay (Qualitative- 100 ng/mL Cutoff Only) REF 100248
Impostazioni DxC 500 AU / 500i Urine Nome calibratore: DRI Oxycodone Calibrator REF 100250

ID reagente 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 ▼
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
REACTION SLOPE	+										
MEASURING POINT	Point 1: First	17		Last	21		LIH INFLUENCE CHECK	☐ Perform LIH check			
	Point 2: First			Last			Lipemia	+	▼		
							Icterus	+	▼		
							Hemolysis	+	▼		
Linearity Limit			%								
Lag Time Check	☐ Perform Lag Time Check										

Nome reagente: DRI Oxycodone Assay (Qualitative- 100 ng/mL Cutoff Only) REF 100248
Nome del calibratore per impostazione urine DxC 500 AU / 500i DRI Oxycodone Calibrator
REF 100250, Continua

ID reagente 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

Add

Positive Cutoff

☐ SLOPE CHECK

Slope Check

Number of Levels

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-2	100	-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	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

Nome reagente: DRI Oxycodone Assay (Qualitative- 300 ng/mL Cutoff Only) REF 100248
Impostazioni DxC 500 AU / 500i Urine Nome calibratore: DRI Oxycodone Calibrator REF
100251

ID reagente 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						Last: Low	-2.0000	High	3.0000
	R1-1	87 μL	Dilution	0 μL	ANALYTICAL MEASURING RANGE	Low	0.00	High	1000.00
	R2-1	87 μL	Dilution	0 μL					
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								

Nome reagente: DRI Oxycodone Assay (Qualitative- 300 ng/mL Cutoff Only) REF 100248
Impostazioni DxC 500 AU / 500i Urine Nome calibratore: DRI Oxycodone Calibrator REF
100251, *Continua*

ID reagente 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

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

	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	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

Nome reagente: DRI Oxycodone Assay (Semi-Quantitative- 100 ng/mL Cutoff Only)

ID reagente 557

REF 100248 Impostazioni DxC 500 AU / 500i Urine

Nome calibratore: DRI Oxycodone Calibrators 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											

Nome reagente: DRI Oxycodone Assay (Semi-Quantitative- 100 ng/mL Cutoff Only)

ID reagente 557

REF 100248 Impostazioni DxC 500 AU / 500i Urine

Nome calibratore DRI Oxycodone Calibrators REF 1664, 100250, 100251, 100252, 100253, *Continua*

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
Add **OXY**

Positive Cutoff **100**

☒ SLOPE CHECK

Number of Levels **5**

Slope Check **+**

STABILITY AND INTERVAL

Reagent Blank Stability Days Hours

Calibration Stability Days Hours

Interval **Bottle**

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**

Nome reagente: DRI Oxycodone Assay (Semi-Quantitative- 300 ng/mL Cutoff Only)

ID reagente 557

REF 100248 Impostazioni DxC 500 AU / 500i Urine

Nome calibratore: DRI Oxycodone Calibrators 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	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										

Nome reagente: DRI Oxycodone Assay (Semi-Quantitative- 300 ng/mL Cutoff Only)

ID reagente 557

REF 100248 Impostazioni DxC 500 AU / 500i Urine

Nome calibratore: DRI Oxycodone Calibrators REF 1664, 100250, 100251, 100252, 100253, *Continua*

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

Limit Points

Limit 1	0
Limit 2	27

Check Pattern

Pattern **Pattern 1**

☐ Logic Check 2

Check Points

Point 1	0
Interval	1

Limit Points

Limit 1	0
Limit 2	27

☐ Logic Check 3

Check Points

Point 1	0
Interval	1

Limit Points

Limit 1	0
Limit 2	27

Decision Values

Value 1	0
Value 2	0

Informazioni supplementari

Importante

Poiché Beckman Coulter non produce il reagente né esegue controlli di qualità o altre analisi sui singoli lotti, Beckman Coulter non può essere responsabile della qualità dei dati ottenuti in base alle prestazioni del reagente, di eventuali variazioni tra i lotti di reagente o delle modifiche al protocollo da parte del produttore.

Danni dovuti al trasporto

Se il prodotto risulta danneggiato alla consegna, contattare il centro di supporto tecnico Beckman Coulter.

© 2024 Thermo Fisher Scientific Inc. Tutti i diritti riservati. Tutti i marchi sono di proprietà di Thermo Fisher Scientific e delle sue consociate salvo diversamente specificato. I sistemi della serie AU sono marchi di Beckman Coulter.



 B·R·A·H·M·S GmbH, Neuendorfstrasse 25, 16761, Hennigsdorf, Germania

Fine