FORM 3

UNITED STATES SECURITIES AND EXCHANGE COMMISSION

Washington, D.C. 20549

INITIAL STATEMENT OF BENEFICIAL OWNERSHIP OF SECURITIES

OMB APPROVAL							
OMB Number: 3235-0104							
Estimated average burden							
hours per response	e: 0.5						

Filed pursuant to Section 16(a) of the Securities Exchange Act of 1934 or Section 30(h) of the Investment Company Act of 1940

1. Name and Ad Connolly J	dress of Reportin	R (N	Date of Event equiring Statem Month/Day/Year 2/06/2017	nent (ame and Ticker or Trac rix Corp [QTRX				
	(First) (Middle) CAPITAL VENTURE			Relationship of Reporting Person(s) to Issuer (Check all applicable) X Director 10% Owner			(5. If Amendment, Date of Original Filed (Month/Day/Year)		
INVESTORS, LLC 200 CLARENDON STREET				Officer (give title below)	Other (spe below)	· ' '	6. Individual or Joint/Group Filing (Check Applicable Line) X Form filed by One Reporting Person			
(Street) BOSTON	MA	02116								y More than One
(City)	(State)	(Zip)								
		T	able I - Non	-Derivati	ve Secu	rities Beneficiall	y Owned			
1. Title of Secui	rity (Instr. 4)	Т	able I - Non	2.	Amount o	rities Beneficiall of Securities Owned (Instr. 4)	3. Ownersh Form: Direct or Indirect (Instr. 5)	cṫ (D) (II	Nature of Indirect	Beneficial Ownership
1. Title of Secu	rity (Instr. 4)		Table II - D	2. Be	Amount o	of Securities	3. Ownersh Form: Direct or Indirect (Instr. 5)	ct (D) (II (I)		Beneficial Ownership
	rity (Instr. 4)	(e.g	Table II - D	erivative S, warrar	Amount of eneficially Security Security 3. Title a	of Securities Owned (Instr. 4)	3. Ownersh Form: Direct or Indirect (Instr. 5) Owned securitie	ct (D) (II (I)	5. Ownership	6. Nature of Indirect Beneficial Ownership (Instr. 5)

Explanation of Responses:

Remarks:

No securities are beneficially owned.

/s/ John M. Connolly 12/06/2017

** Signature of Reporting Person Date

Reminder: Report on a separate line for each class of securities beneficially owned directly or indirectly.

Note: File three copies of this Form, one of which must be manually signed. If space is insufficient, see Instruction 6 for procedure.

Persons who respond to the collection of information contained in this form are not required to respond unless the form displays a currently valid OMB Number.

^{*} If the form is filed by more than one reporting person, see Instruction 5 (b)(v).

^{**} Intentional misstatements or omissions of facts constitute Federal Criminal Violations See 18 U.S.C. 1001 and 15 U.S.C. 78ff(a).